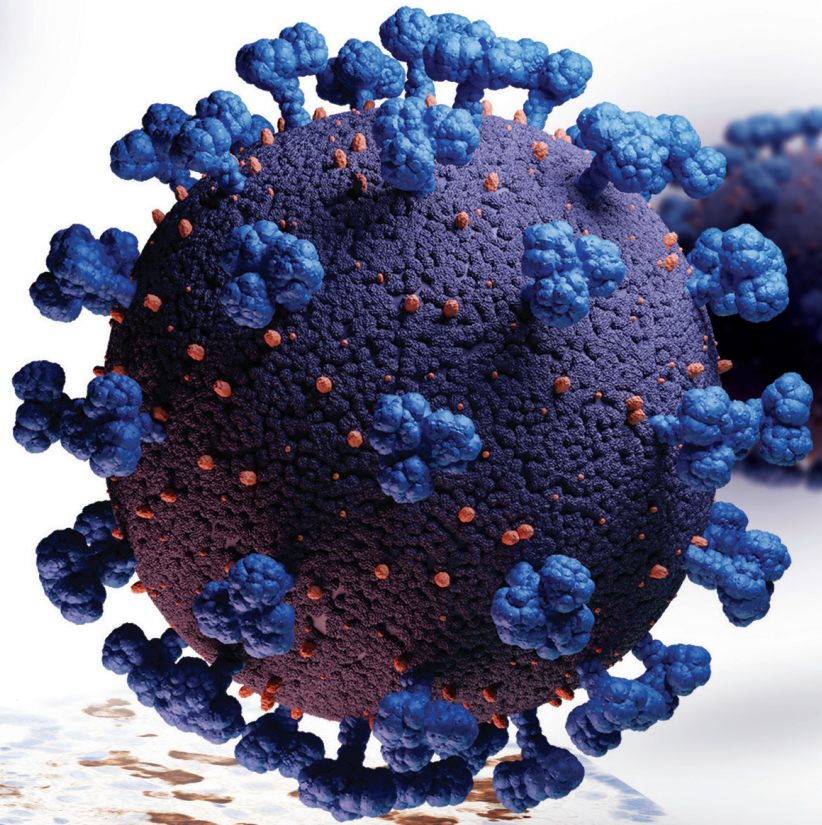


Powered by
RNAscope™
RNA *In Situ* Hybridization

RNAscope ISH Probes for Anatomic Pathology



biotechne®



Superior Sensitivity & Specificity

From our wide portfolio of ready-to-use probes, ACD provides a broad menu of high-quality RNAscope™ ISH probes for many pathological RNA markers. Using our proprietary ACD RNAscope Probe Design Pipeline, we provide double “Z” oligo probes designed to hybridize to your specific RNA target of interest. The RNAscope RNA ISH probes provide labs with the opportunity to add new markers to complement and enhance their traditional IHC applications.

Why RNAscope RNA ISH?

- Unparalleled sensitivity and specificity—
Detect any marker of interest to complement your IHC or DNA ISH, with consistently reliable results.
- Surpass traditional IHC challenges—
Easy validation, omit issues with high background, low signal or suboptimal antibodies.
- Choose your platform—RNAscope RNA ISH probes are available and ready-to-use in both manual and automated formats. RNAscope probes have been developed for full automation on the Leica BOND (LS) and Ventana Discovery (VS) platforms.

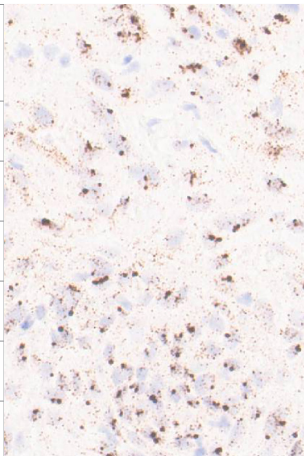
Target Gene	Probe Name	Probe Type	Page
UBC Ubiquitin C	Hs-UBC	Positive control	6
PPIB Peptidylprolyl isomerase B	Hs-PPIB	Positive control	6
POLR2A RNA polymerase II subunit A	Hs-POLR2A	Positive control	6
dapB <i>Bacillus subtilis</i> 4-hydroxy-tetrahydrodipicolinate reductase	DapB	Negative control	7
HBZ T-lymphotropic virus 1 bZIP factor	V-HTVL-HBZ	Target probe	7
E6/E7 <i>Human papillomavirus (HPV) type 16</i>	HPV16	Target probe	7
E6/E7 HPV18	HPV18	Target probe	7
E6/E7 HPV31	HPV31	Target probe	8
E6/E7 HPV33	HPV33	Target probe	8
E6/E7 HPV45	HPV45	Target probe	8
E6/E7 HPV16/18	HPV16/18	Target probe	8
E6/E7 HPV High Risk 18 subtypes	HPV-HR18	Target probe	9
E6/E7 HPV Low Risk 6 subtypes	HPV LR6	Target probe	9
16S rRNA <i>Mycobacterium tuberculosis</i> H37Rv	B-MTB-NTM-16SrRNA-O1	Target probe	9
CDKN2A Cyclin-dependent kinase inhibitor 2A	Hs-CDKN2A	Target probe	9
ESR1 Estrogen receptor 1	Hs-ESR1	Target probe	10
CXCL13 C-X-C motif chemokine ligand 13	Hs-CXCL13	Target probe	10
MDM2 oncogene E3 ubiquitin protein ligase	Hs-MDM2	Target probe	10
ALK Alk receptor tyrosine kinase	Hs-ALK	Target probe	10
IGKC Immunoglobulin kappa locus constant region	Hs-IGK	Target probe	11
IGLC1 Immunoglobulin lambda locus constant region	Hs-IGL	Target probe	11
IGLL5 Immunoglobulin lambda-like polypeptide 5	Hs-IGLL5	Target probe	11
23S rRNA <i>Mycobacterium tuberculosis</i> H37Rv	B-MTB-23SrRNA-1-C1	Target probe	11
CSF1 Colony stimulating factor 1	Hs-CSF1	Target probe	12
PRAME Preferentially expressed antigen in melanoma	Hs-PRAME	Target probe	12
NAPSA Napsin A aspartic peptidase	Hs-NAPSA	Target probe	12
PDGFB platelet-derived growth factor beta transcript variant 2	Hs-PDGFB	Target probe	12
PDGFB platelet-derived growth factor beta transcript variant 1	Hs-PDGFB-No-XMm	Target probe	13
TNFSF11 TNF (ligand) superfamily member 11	Hs-TNFSF11	Target probe	13
TNFRSF11A TNF receptor superfamily member 11a	Hs-TNFRSF11A	Target probe	13
S Spike glycoprotein SARS-CoV-2 virus	RNAscope ISH Probe SARS-CoV-2	Target probe	13
S Spike glycoprotein SARS-CoV-2 virus	V-nCoV2019-S	Target probe	13
MYB MYB proto-oncogene	Hs-MYB	Target probe	14
NOS2 Nitric oxide synthase 2	Hs-NOS2	Target probe	14
AFP Alpha fetoprotein	Hs-AFP	Target probe	14

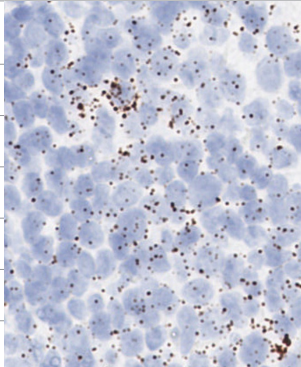
Target Gene	Probe Name	Probe Type	Page
HAVCR1 Hepatitis A virus (<i>HAV</i>) cellular receptor 1	Hs-HAVCR1	Target probe	14
FCRL4 Fc receptor-like 4	Hs-FCRL4	Target probe	15
KRT17 Keratin 17	Hs-KRT17	Target probe	15
PDCD1LG2 Programmed cell death 1 ligand 2	Hs-PDCD1LG2	Target probe	15
FGF23 Fibroblast growth factor 23	Hs-FGF23	Target probe	15
ALB Albumin	Hs-ALB	Target probe	16
TERT Telomerase reverse transcriptase	Hs-TERT	Target probe	16
18S rRNA <i>Candida albican</i>	F-C.albicans-18SrRNA	Target probe	16
EBER-1 Epstein-Barr virus (<i>EBV</i>)	V-EBV-EBER1	Target probe	16
RNA2.7 Human cytomegalovirus (<i>CMV</i>)	V-CMV	Target probe	17
23S rRNA <i>Bartonella henselae</i> strain Houston-1	B-B.henselae-23SrRNA	Target probe	17
16S rRNA <i>Helicobacter pylori</i> strain ATCC 4350	B-H.pylori-16S	Target probe	17
23S rRNA <i>Treponema pallidum</i> subsp. pallidum strain SS14	B-T.pallidum-23SrRNA	Target probe	17
NP Nucleoprotein Respiratory syncytial virus	V-RSV-NP	Target probe	18
LargeT BK polyomavirus (<i>BKPyV</i>)	V-BKPyV-largeT	Target probe	18
LargeAg Merkel cell polyomavirus isolate R17b (<i>MCPyV</i>)	V-MCPyV-LT-ST-Ag-O1	Target probe	18

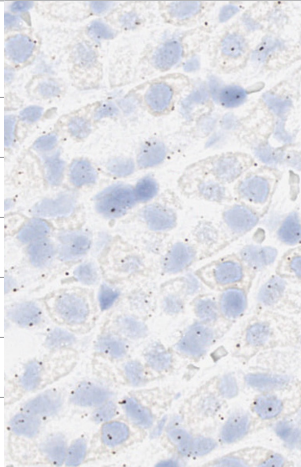
The species is typically identified by one or two letter designations such as "B" for Bacteria, "F" for Fungi, "V" for Virus, and "Hs" for Homo Sapiens.

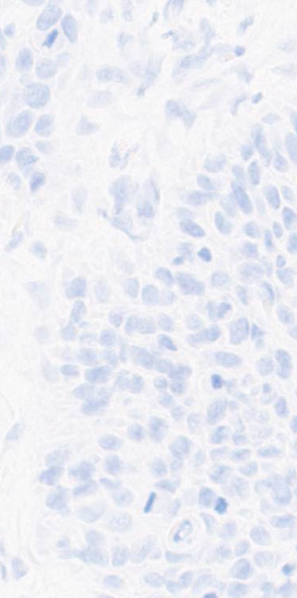
Portfolio of ready-to-use probes

Intuitive protocol, implemented seamlessly into your workflow
 - The RNAscope RNA ISH protocol is easy to follow, similar in many steps to IHC and can be transferred to your established Leica BOND (LS) or Ventana Discovery (VS) staining platform with ease.

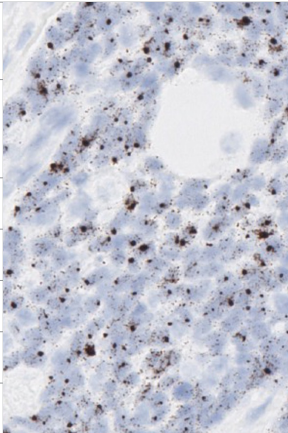
Hs-UBC (Positive Control)		
Catalog No.	310041	RNAscope Positive Control Probe - Hs-UBC
	312028	RNAscope 2.5 LS Positive Control Probe - Hs-UBC
	312029	RNAscope 2.5 VS Positive Control Probe - Hs-UBC
Probe Description	Positive control probe for medium high level expressing targets. For Research Use Only.† UBC positive control probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	Homo sapiens ubiquitin C (UBC) mRNA	
Species	Homo sapiens	
Target Gene	UBC	
Entrez Gene ID	7316	
Accession#	NM_021009	
Probe Region	342 - 1503	
ZZ Pair No.	20	

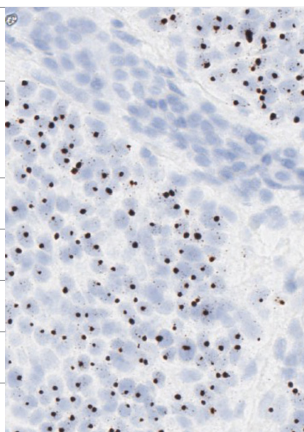
Hs-PPIB (Positive Control)		
Catalog No.	313901	RNAscope Positive Control Probe - Hs-PPIB
	313908	RNAscope 2.5 LS Positive Control Probe_Hs-PPIB
	313909	RNAscope 2.5 VS Positive Control Probe_Hs-PPIB
Probe Description	Positive control probe for medium level expressing targets. For Research Use Only.† PPIB positive control probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	Homo sapiens PPIB mRNA	
Species	Homo sapiens	
Target Gene	PPIB	
Entrez Gene ID	5479	
Accession#	NM_000942.4	
Probe Region	138-989	
ZZ Pair No.	16	

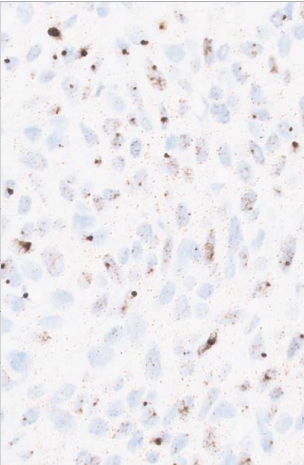
Hs-POLR2A (Positive Control)		
Catalog No.	310451	RNAscope Positive Control Probe - Hs-POLR2A
	310458	RNAscope 2.5 LS Probe - Hs-POLR2A
	310459	RNAscope 2.5 VS Positive Control Probe - Hs-POLR2A
Probe Description	Positive control probe for low level expressing targets. For Research Use Only.†	
Specificity	Homo sapiens POLR2A, 220kDa mRNA	
Species	Homo sapiens	
Target Gene	POLR2A	
Entrez Gene ID	5430	
Accession#	NM_000937.4	
Probe Region	2514-3433	
ZZ Pair No.	20	

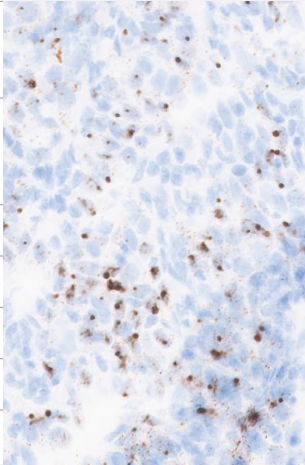
dapB (Negative Control)		
Catalog No.	310043	RNAscope Negative Control Probe - DapB
	312038	RNAscope 2.5 LS Negative Control Probe_dapB
	312039	RNAscope 2.5 VS Negative Control Probe_dapB
Probe Description	Negative control probe. For Research Use Only.† dapB negative control probe is also available as an Analyte Specific Reagent (ASRs) via Leica Biosystems.	
Specificity	<i>Bacillus subtilis</i> subsp. subtilis str. SMY, 4-hydroxy-tetrahydrodipicolinate reductase (dapB) mRNA transcripts.	
Species	<i>Bacillus subtilis</i>	
Target Gene	dapB	
Entrez Gene ID	939024	
Accession#	EF191515	
Probe Region	414-862	
ZZ Pair No.	10	

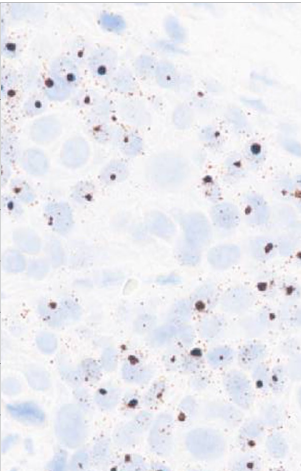
V-HTLV-HBZ		
Catalog No.	432901	RNAscope Probe- V-HT-VL-HBZ
	432908	RNAscope 2.5 LS Probe- V-HT-VL-HBZ
	432909	RNAscope 2.5 VS Probe - V-HT-VL-HBZ
Probe Description	Target probe for Human T-lymphotropic virus 1 bZIP factor (HBZ) mRNA. For Research Use Only.†	
Specificity	HTLV HBZ mRNA transcript	
Species	Human T-lymphotropic virus 1	
Target Gene	HBZ	
Entrez Gene ID	N/A	
Accession#	DQ273132.1	
Probe Region	173 - 1363	
ZZ Pair No.	20	

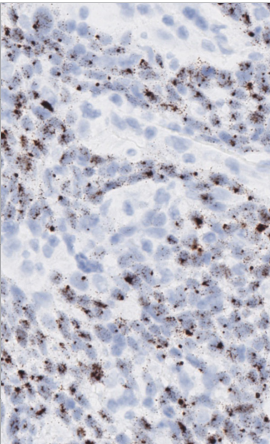
HPV 16		
Catalog No.	311521	RNAscope Probe - HPV16
	311528	RNAscope 2.5 LS Probe - HPV16
	311529	RNAscope 2.5 VS Probe - HPV16
Probe Description	Target probe for human papillomavirus type 16 E6/E7 mRNA. For Research Use Only.† HPV16 target probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	HPV type 16 E6/E7 mRNA	
Species	Human papillomavirus type 16	
Target Gene	E6, E7	
Entrez Gene ID	N/A	
Accession#	KF880690.1	
Probe Region	84-753	
ZZ Pair No.	10	

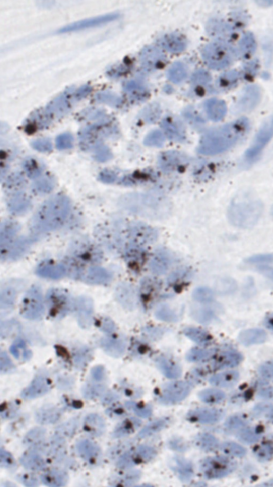
HPV18		
Catalog No.	311531	RNAscope Probe - HPV18
	311538	RNAscope 2.5 LS Probe - HPV18
	311539	RNAscope 2.5 VS Probe - HPV18
Probe Description	Target probe for human papillomavirus type 18 E6/E7 mRNA. For Research Use Only.† HPV18 target probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	HPV type 18 E6/E7 mRNA	
Species	Human papillomavirus type 18	
Target Gene	E6, E7	
Entrez Gene ID	N/A	
Accession#	AY262282.1	
Probe Region	142-838	
ZZ Pair No.	10	

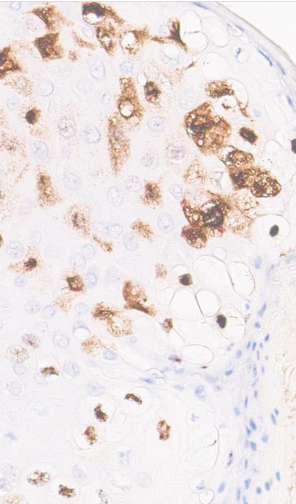
HPV31		
Catalog No.	311551	RNAscope Probe - HPV31
	311558	RNAscope 2.5 LS Probe - HPV31
	311559	RNAscope 2.5 VS Probe - HPV31
Probe Description	Target probe for human papillomavirus type 31 E6/E7 mRNA. For Research Use Only.† HPV31 target probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	HPV type 31 E6/E7 mRNA	
Species	Human papillomavirus type 31	
Target Gene	E6, E7	
Entrez Gene ID	N/A	
Accession#	J04353.1	
Probe Region	110-755	
ZZ Pair No.	10	

HPV33		
Catalog No.	311561	RNAscope Probe - HPV33
	311568	RNAscope 2.5 LS Probe - HPV33
	311569	RNAscope 2.5 VS Probe - HPV33
Probe Description	Target probe for human papillomavirus type 33 E6/E7 mRNA. For Research Use Only.† HPV33 target probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	HPV type 33 E6/E7 mRNA	
Species	Human papillomavirus type 33	
Target Gene	E6, E7	
Entrez Gene ID	N/A	
Accession#	M12732.1	
Probe Region	114-840	
ZZ Pair No.	10	

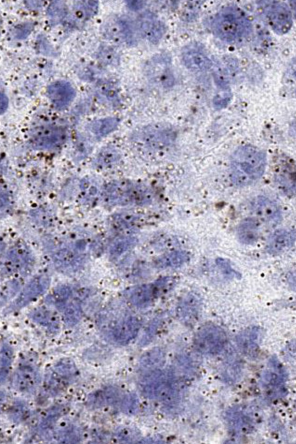
HPV45		
Catalog No.	311591	RNAscope Probe - HPV45
	311598	RNAscope 2.5 LS Probe - HPV45
	311599	RNAscope 2.5 VS Probe - HPV45
Probe Description	Target probe for human papillomavirus type 45 E6/E7 mRNA. For Research Use Only.†	
Specificity	HPV type 45 E6/E7 mRNA	
Species	Human papillomavirus type 45	
Target Gene	E6, E7	
Entrez Gene ID	N/A	
Accession#	X74479.1	
Probe Region	140-838	
ZZ Pair No.	10	

HPV16/18		
Catalog No.	311121	RNAscope 2.5 Probe - HPV16/18
	311128	RNAscope 2.5 LS Probe - HPV16/18
	311129	RNAscope 2.5 VS Probe - HPV16/18
Probe Description	Pooled target probes for HPV high-risk type 16 and 18 E6/E7 mRNA. For Research Use Only.†	
Specificity	HPV high-risk type 16 and 18 E6/E7 mRNA	
Species	Human papillomavirus type 16 and 18	
Target Gene	E6, E7	
Entrez Gene ID	NA	
Accession#	NA	
Probe Region	NA	
ZZ Pair No.	NA	

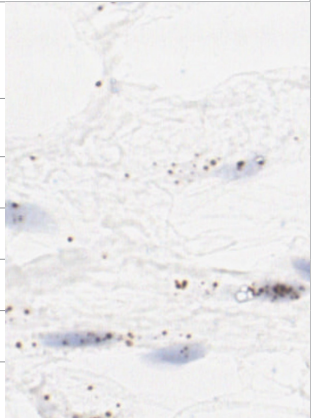
HPV HR18		
Catalog No.	312591	RNAscope Probe - HPV-HR18
	312598	RNAscope 2.5 LS Probe - HPV-HR18
	312599	RNAscope 2.5 VS Probe - HPV-HR18
Probe Description	Pooled target probes for HPV high-risk type 16, 18, 26, 31, 33, 35, 39, 45, 51, 52, 53, 56, 58, 59, 66, 68, 73 and 82, E6/E7 mRNA. For Research Use Only.†	
Specificity	HPV high-risk type 16, 18, 26, 31, 33, 35, 39, 45, 51, 52, 53, 56, 58, 59, 66, 68, 73 and 82, E6/E7 mRNA.	
Species	Human papillomavirus	
Target Gene	E6, E7	
Entrez Gene ID	NA	
Accession#	NA	
Probe Region	NA	
ZZ Pair No.	NA	

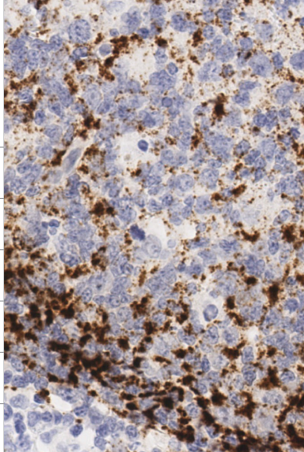
HPV LR6		
Catalog No.	407601	RNAscope Probe - HPV LR6
	407608	RNAscope 2.5 LS Probe - HPV LR6
	407609	RNAscope 2.5 VS Probe - HPV LR6
Probe Description	Pooled target probes for HPV low-risk type 6, 11, 40, 42, 43, 44 E6/E7 mRNA. For Research Use Only.†	
Specificity	HPV low-risk type 6, 11, 40, 42, 43, 44 E6/E7 mRNA.	
Species	Human papillomavirus	
Target Gene	E6, E7	
Entrez Gene ID	N/A	
Accession#	N/A	
Probe Region	N/A	
ZZ Pair No.	N/A	

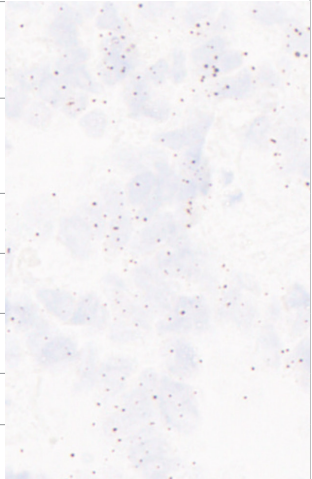
B-MTB-NTM-16SrRNA-O1		
Catalog No.	496191	RNAscope Probe - B-MTB-NTM-16SrRNA-O1
	496198	RNAscope 2.5 LS Probe-B-MTB-NTM-16SrRNA-O1
	496199	RNAscope 2.5 VS Probe-B-MTB-NTM-16SrRNA-O1
Probe Description	Target probe for <i>Mycobacterium tuberculosis</i> H37Rv 16S ribosomal RNA. For Research Use Only.†	
Specificity	MTB 16S rRNA	
Species	<i>Mycobacterium tuberculosis</i>	
Target Gene	16S rRNA	
Entrez Gene ID	N/A	
Accession#	N/A	
Probe Region	1-40	
ZZ Pair No.	1	

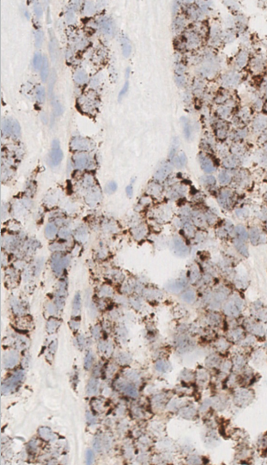
Hs-CDKN2A		
Catalog No.	310181	RNAscope Probe - Hs-CDKN2A
	310188	RNAscope LS 2.5 Probe - Hs-CDKN2A
	310189	RNAscope 2.5 VS Probe - Hs-CDKN2A
Probe Description	Target probe for Homo sapiens cyclin-dependent kinase inhibitor 2A (CDKN2A) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens CDKN2A transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	CDKN2A	
Entrez Gene ID	1029	
Accession#	NM_000077.4	
Probe Region	95 - 1206	
ZZ Pair No.	17	

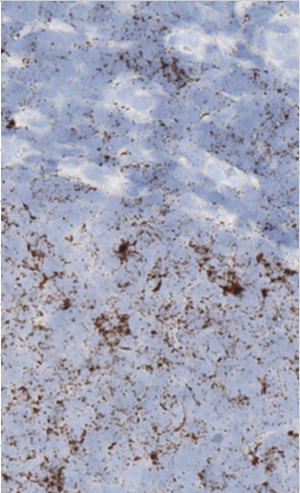
Hs-ESR1		
Catalog No.	310301	RNAscope Probe - Hs-ESR1
	310308	RNAscope LS 2.5 Probe - Hs-ESR1
	310309	RNAscope 2.5 VS Probe - Hs-ESR1
Probe Description	Target probe for Homo sapiens estrogen receptor 1 (ESR1) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens ESR1 transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	ESR1	
Entrez Gene ID	2099	
Accession#	NM_000125.3	
Probe Region	1251-2376	
ZZ Pair No.	20	

Hs-MDM2		
Catalog No.	311781	RNAscope Probe - Hs-MDM2
	311788	RNAscope 2.5 LS Probe - Hs-MDM2
	311789	RNAscope 2.5 VS Probe - Hs-MDM2
Probe Description	Target probe for Homo sapiens MDM2 oncogene, E3 ubiquitin protein ligase (MDM2) mRNA. For Research Use Only.†	
Specificity	Homo sapiens MDM2 mRNA	
Species	Homo sapiens	
Target Gene	MDM2	
Entrez Gene ID	4193	
Accession#	NM_002392.4	
Probe Region	289-1511	
ZZ Pair No.	20	

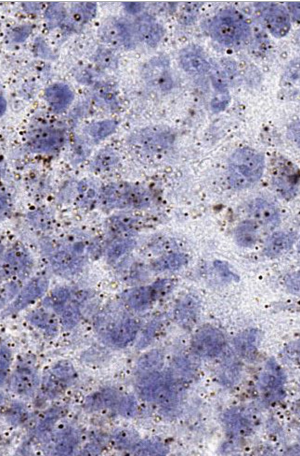
Hs-CXCL13		
Catalog No.	311321	RNAscope Probe- Hs-CXCL13
	311328	RNAscope LS 2.5 Probe- Hs-CXCL13
	311329	RNAscope VS 2.5 Probe- Hs-CXCL13
Probe Description	Target probe for Homo sapiens C-X-C motif chemokine ligand 13 (CXCL13) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens CXCL13 transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	CXCL13	
Entrez Gene ID	10563	
Accession#	NM_006419.2	
Probe Region	30 - 1186	
ZZ Pair No.	16	

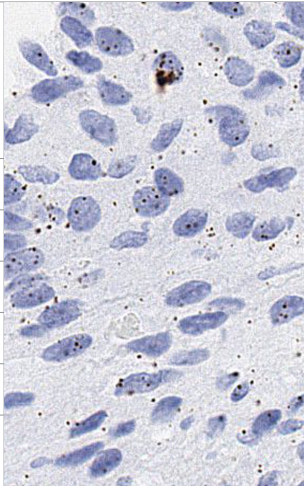
Hs-ALK		
Catalog No.	311841	RNAscope Probe - Hs-ALK
	311848	RNAscope LS 2.5 Probe - Hs-ALK
	311849	RNAscope 2.5 VS Probe - Hs-ALK
Probe Description	Target probe for Homo sapiens ALK receptor tyrosine kinase (ALK) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens ALK mRNA	
Species	Homo sapiens	
Target Gene	ALK	
Entrez Gene ID	238	
Accession#	NM_004304.4	
Probe Region	4749 - 5794	
ZZ Pair No.	20	

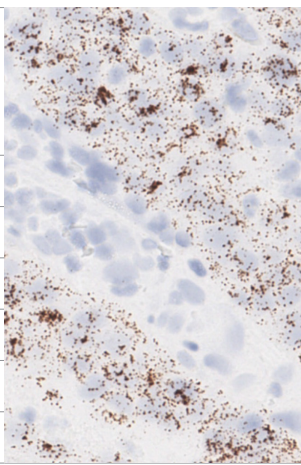
Hs-IGK		
Catalog No.	312551	RNAscope Probe - Hs-IGK
	312558	RNAscope LS 2.5 Probe - Hs-IGK
	312559	RNAscope 2.5 VS Probe - Hs-IGK
Probe Description	Target probe for Human sapiens immunoglobulin kappa locus constant region. For Research Use Only.†	
Specificity	Immunoglobulin kappa locus constant region mRNA transcript	
Species	Homo sapiens	
Target Gene	IGKC	
Entrez Gene ID	3514	
Accession#	N/A	
Probe Region	RC gi 528476665:89086517-89086775	
ZZ Pair No.	6	

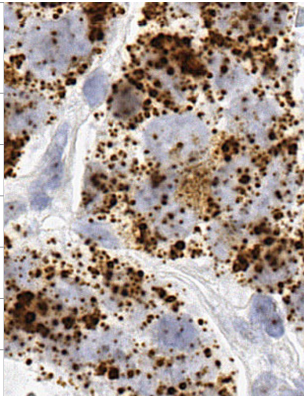
Hs-IGL		
Catalog No.	312561	RNAscope Probe - Hs-IGL
	312568	RNAscope LS 2.5 Probe - Hs-IGL
	312569	RNAscope 2.5 VS Probe - Hs-IGL
Probe Description	Target probe for Human sapiens immunoglobulin lambda locus constant region. For Research Use Only.†	
Specificity	Homo sapiens Immunoglobulin lambda constant region mRNA	
Species	Homo sapiens	
Target Gene	IGLC1	
Entrez Gene ID	3537	
Accession#	N/A	
Probe Region	gi 528476531:23255447-23255726	
ZZ Pair No.	7	

B-MTB-23SrRNA		
Catalog No.	471651	RNAscope Probe - B-MTB-23SrRNA-1-C1
	471658	RNAscope 2.5 LS Probe - B-MTB-23SrRNA-1-C1
	471659	RNAscope 2.5 VS Probe - B-MTB-23SrRNA-1-C1
Probe Description	Target probe for <i>Mycobacterium tuberculosis</i> H37Rv 23S ribosomal RNA. For Research Use Only.†	
Specificity	MTB 23S rRNA	
Species	<i>Mycobacterium tuberculosis</i> H37Rv	
Target Gene	23S rRNA	
Entrez Gene ID	N/A	
Accession#	AL123456	
Probe Region	1474379-1474455	
ZZ Pair No.	1	

Hs-IgLL5		
Catalog No.	410271	RNAscope Probe - Hs-IgLL5
	410278	RNAscope LS 2.5 Probe - Hs-IgLL5
	410279	RNAscope 2.5 VS Probe - Hs-IgLL5
Probe Description	Target probe for Homo sapiens immunoglobulin lambda-like polypeptide 5 (IgLL5) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens IgLL5 transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	IgLL5	
Entrez Gene ID	100423062	
Accession#	NM_001178126.1	
Probe Region	37 - 448	
ZZ Pair No.	6	

Hs-CSF1		
Catalog No.	313001	RNAscope Probe - Hs-CSF1
	313008	RNAscope LS 2.5 Probe - Hs-CSF1
	313009	RNAscope VS 2.5 Probe - Hs-CSF1
Probe Description	Target probe for Homo sapiens colony stimulating factor 1 (macrophage) (CSF1) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens CSF1 transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	CSF1	
Entrez Gene ID	1435	
Accession#	NM_000757.5	
Probe Region	Feb-68	
ZZ Pair No.	30	

Hs-PRAME		
Catalog No.	401851	RNAscope Probe - Hs-PRAME
	401858	RNAscope LS 2.5 Probe - Hs-PRAME
	401859	RNAscope 2.5 VS Probe - Hs-PRAME
Probe Description	Target probe for Homo sapiens preferentially expressed antigen in melanoma (PRAME) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens PRAME transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	PRAME	
Entrez Gene ID	23532	
Accession#	NM_006115.3	
Probe Region	760-1811	
ZZ Pair No.	20	

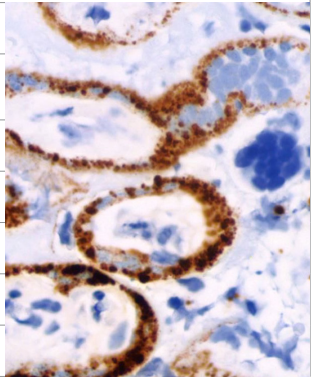
Hs-NAPSA		
Catalog No.	405821	RNAscope Probe - Hs-NAPSA
	405828	RNAscope LS 2.5 Probe - Hs-NAPSA
	405829	RNAscope 2.5 VS Probe - Hs-NAPSA
Probe Description	Target probe for Homo sapiens napsin A aspartic peptidase (NAPSA) mRNA. For Research Use Only.† Napsin A target probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	Homo sapiens NAPSA mRNA	
Species	Homo sapiens	
Target Gene	NAPSA	
Entrez Gene ID	9476	
Accession#	NM_004851.1	
Probe Region	2-895	
ZZ Pair No.	20	

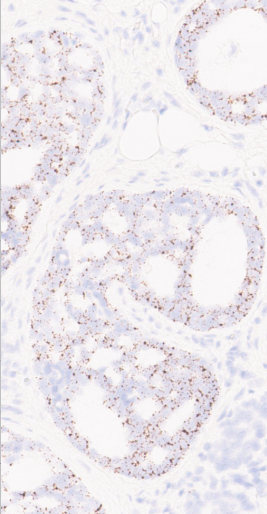
Hs-PDGFB		
Catalog No.	406701	RNAscope Probe - Hs-PDGFB
	406708	RNAscope LS 2.5 Probe- Hs-PDGFB
	406709	RNAscope VS 2.5 Probe- Hs-PDGFB
Probe Description	Target probe for Homo sapiens platelet-derived growth factor beta polypeptide (PDGFB) transcript variant 2 mRNA. For Research Use Only.†	
Specificity	PDGFB transcript variant 2 mRNA	
Species	Homo sapiens	
Target Gene	PDGFB	
Entrez Gene ID	5155	
Accession#	NM_033016.2	
Probe Region	665 - 2037	
ZZ Pair No.	21	

Hs-PDGFB-No-XMm		
Catalog No.	441571	RNAscope Probe - Hs-PDGFB-No-XMm
	441578	RNAscope LS 2.5 Probe - Hs-PDGFB-No-XMm
	441579	RNAscope 2.5 VS Probe - Hs-PDGFB-No-XMm
Probe Description	Target probe for Homo sapiens platelet-derived growth factor beta polypeptide (PDGFB) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens PDGFB transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	PDGFB	
Entrez Gene ID	5155	
Accession#	NM_002608.2	
Probe Region	1995-6047	
ZZ Pair No.	20	

Hs-TNFSF11		
Catalog No.	412271	RNAscope Probe- Hs-TNFSF11
	412278	RNAscope LS 2.5 Probe- Hs-TNFSF11
	412279	RNAscope VS 2.5 Probe- Hs-TNFSF11
Probe Description	Target probe for Homo sapiens tumor necrosis factor (ligand) superfamily member 11 (TNFSF11) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens TNFSF11 transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	TNFSF11	
Entrez Gene ID	8600	
Accession#	NM_003701.3	
Probe Region	366-1507	
ZZ Pair No.	20	

Hs-TNFRSF11A		
Catalog No.	412281	RNAscope Probe- Hs-TNFRSF11A
	412288	RNAscope 2.5 LS Probe- Hs-TNFRSF11A
	412289	RNAscope 2.5 VS Probe- Hs-TNFRSF11A
Probe Description	Target probe for Homo sapiens tumor necrosis factor receptor superfamily member 11a NFKB activator (TNFRSF11A) transcript variant 4 mRNA. For Research Use Only.†	
Specificity	Homo sapiens TNF receptor superfamily member 11a (TNFRSF11A) transcript variant 4 mRNA	
Species	Homo sapiens	
Target Gene	TNFRSF11A	
Entrez Gene ID	8792	
Accession#	NM_001270951.1	
Probe Region	1928 - 2981	
ZZ Pair No.	20	

SARS-CoV-2		
Catalog No.	848561	RNAscope Probe - V-nCoV2019-S
	848568	RNAscope 2.5 LS Probe - V-nCoV2019-S
	848569	RNAscope 2.5 VS Probe - V-nCoV2019-S
	200440*	RNAscope ISH Probe SARS-CoV-2 (ASR)
Probe Description	Target probe for severe acute respiratory syndrome coronavirus 2 isolate Wuhan-Hu-1 spike glycoprotein (S) mRNA. For Research Use Only.† The RNAscope SARS-CoV-2 probe is also available as an Analyte Specific Reagent (ASR). <i>* Analyte Specific Reagent. Analytical and performance characteristics are not established.</i>	
Specificity	S gene mRNA	
Species	SARS-CoV-2 virus	
Target Gene	S	
Entrez Gene ID	N/A	
Accession#	NC_045512.2	
Probe Region	21631 - 23303	
ZZ Pair No.	20	

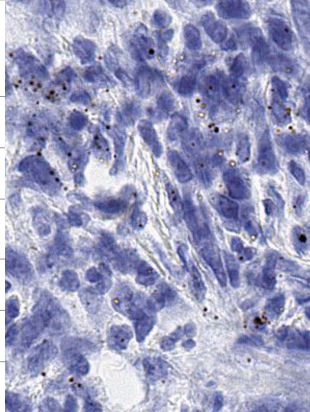
Hs-MYB		
Catalog No.	419701	RNAscope Probe - Hs-MYB
	419708	RNAscope LS 2.5 Probe - Hs-MYB
	419709	RNAscope 2.5 VS Probe - Hs-MYB
Probe Description	Target probe for Homo sapiens v-myb avian myeloblastosis viral oncogene homolog (MYB) transcript variant 7 mRNA. For Research Use Only.†	
Specificity	Homo sapiens MYB proto-oncogene, transcription factor (MYB), transcript variant 7 mRNA	
Species	Homo sapiens	
Target Gene	MYB	
Entrez Gene ID	4602	
Accession#	NM_001161659.1t	
Probe Region	204-1289	
ZZ Pair No.	20	

Hs-NOS2		
Catalog No.	424991	RNAscope Probe - Hs-NOS2
	424998	RNAscope LS 2.5 Probe - Hs-NOS2
	424999	RNAscope 2.5 VS Probe - Hs-NOS2
Probe Description	Target probe for Homo sapiens nitric oxide synthase 2 (NOS2) mRNA. For Research Use Only.†	
Specificity	Homo sapiens NOS2 mRNA	
Species	Homo sapiens	
Target Gene	NOS2	
Entrez Gene ID	4843	
Accession#	NM_000625.4	
Probe Region	27-1129	
ZZ Pair No.	20	

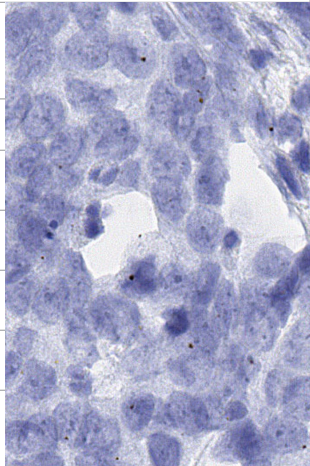
Hs-AFP		
Catalog No.	427411	RNAscope Probe - Hs-AFP
	427418	RNAscope LS 2.5 Probe - Hs-AFP
	427419	RNAscope 2.5 VS Probe - Hs-AFP
Probe Description	Target probe for Homo sapiens alpha fetoprotein (AFP), transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens AFP transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	AFP	
Entrez Gene ID	174	
Accession#	NM_001134.2	
Probe Region	2 - 1078	
ZZ Pair No.	20	

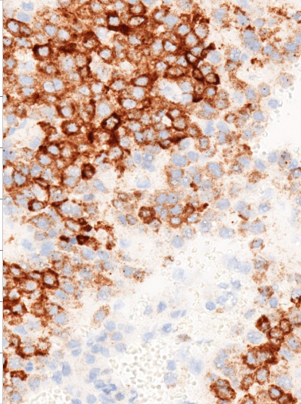
Hs-HAVCR1		
Catalog No.	452141	RNAscope Probe - Hs-HAVCR1
	452148	RNAscope LS 2.5 Probe - Hs-HAVCR1
	452149	RNAscope 2.5 VS Probe - Hs-HAVCR1
Probe Description	Target probe for Homo sapiens hepatitis A virus cellular receptor 1 (HAVCR1) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens HAVCR1 transcripts variant 1 mRNA	
Species	Homo sapiens	
Target Gene	HAVCR1	
Entrez Gene ID	26762	
Accession#	NM_012206.3	
Probe Region	2-1687	
ZZ Pair No.	10	

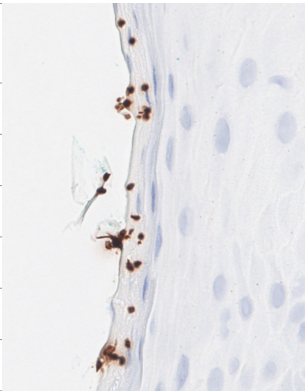
Hs-FCRL4		
Catalog No.	433301	RNAscope Probe - Hs-FCRL4
	433308	RNAscope LS 2.5 Probe - Hs-FCRL4
	433309	RNAscope 2.5 VS Probe - Hs-FCRL4
Probe Description	Target probe for Homo sapiens Fc receptor-like 4 (FCRL4) mRNA. For Research Use Only.†	
Specificity	Homo sapiens FCRL4 mRNA	
Species	Homo sapiens	
Target Gene	FCRL4	
Entrez Gene ID	83417	
Accession#	NM_031282.2	
Probe Region	236-2400	
ZZ Pair No.	30	

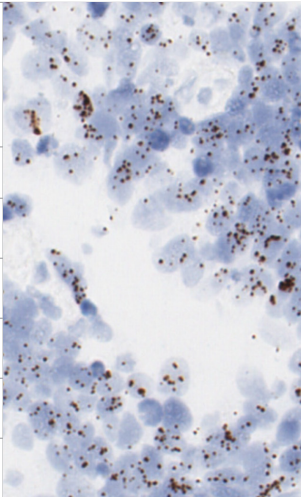
Hs-PDCD1LG2		
Catalog No.	551891	RNAscope Probe- Hs-PDCD1LG2
	551898	RNAscope LS 2.5 Probe- Hs-PDCD1LG2
	551899	RNAscope VS 2.5 Probe- Hs-PDCD1LG2
Probe Description	Target probe for Homo sapiens programmed cell death 1 ligand 2 (PDCD1LG2) mRNA. For Research Use Only.†	
Specificity	Homo sapiens PD-L2 mRNA	
Species	Homo sapiens	
Target Gene	PDCD1LG2	
Entrez Gene ID	80380	
Accession#	NM_025239.3	
Probe Region	222 - 1241	
ZZ Pair No.	20	

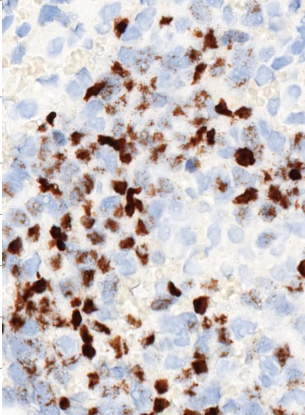
Hs-KRT17		
Catalog No.	463661	RNAscope Probe - Hs-KRT17
	463668	RNAscope 2.5 LS Probe - Hs-KRT17
	463669	RNAscope 2.5 VS Probe - Hs-KRT17
Probe Description	Target probe for Homo sapiens keratin 17 (KRT17) mRNA. For Research Use Only.†	
Specificity	Homo sapiens KRT17 mRNA	
Species	Homo sapiens	
Target Gene	KRT17	
Entrez Gene ID	3872	
Accession#	NM_000422	
Probe Region	302 - 1346	
ZZ Pair No.	19	

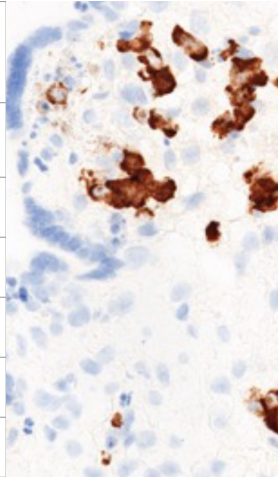
Hs-FGF23		
Catalog No.	557241	RNAscope Probe - Hs-FGF23
	557248	RNAscope LS 2.5 Probe - Hs-FGF23
	557249	RNAscope 2.5 VS Probe - Hs-FGF23
Probe Description	Target probe for Homo sapiens fibroblast growth factor 23 (FGF23) mRNA. For Research Use Only.†	
Specificity	Homo sapiens FGF23 mRNA	
Species	Homo sapiens	
Target Gene	FGF23	
Entrez Gene ID	8074	
Accession#	NM_020638.2	
Probe Region	1275-1988	
ZZ Pair No.	20	

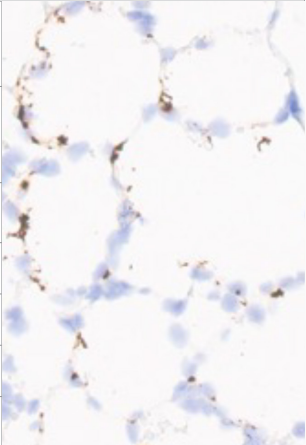
Hs-ALB		
Catalog No.	600941	RNAscope Probe - Hs-ALB
	600948	RNAscope LS 2.5 Probe - Hs-ALB
	600949	RNAscope 2.5 VS Probe - Hs-ALB
Probe Description	Target probe for Homo sapiens albumin (ALB) mRNA. For Research Use Only.† Albumin target probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	Homo sapiens albumin mRNA	
Species	Homo sapiens	
Target Gene	ALB	
Entrez Gene ID	213	
Accession#	NM_000477.5	
Probe Region	164-1183	
ZZ Pair No.	20	

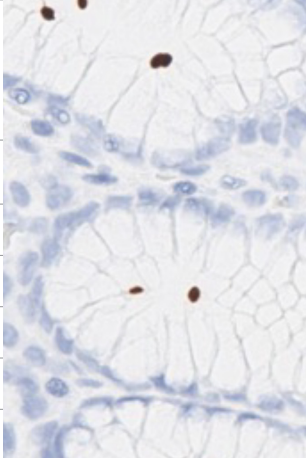
F-C.albicans - 18SrRNA		
Catalog No.	432111	RNAscope Probe - F-C. albicans-18SrRNA
	432118	RNAscope LS 2.5 Probe- F-C. albicans-18SrRNA
	432119	RNAscope 2.5 VS Probe- F-C.albicans-18SrRNA
Probe Description	Target probe for <i>Candida albicans</i> 18S ribosomal RNA. For Research Use Only.†	
Specificity	C. albicans 18S rRNA	
Species	C. albicans	
Target Gene	18S rRNA	
Entrez Gene ID	N/A	
Accession#	E15168.1	
Probe Region	625-1088	
ZZ Pair No.	2	

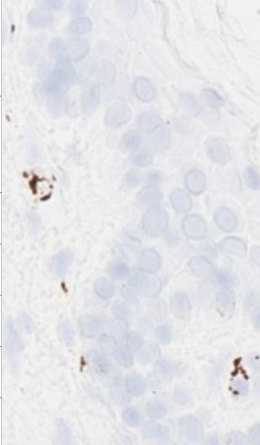
Hs-TERT		
Catalog No.	605511	RNAscope Probe - Hs-TERT
	605518	RNAscope LS 2.5 Probe - Hs-TERT
	605519	RNAscope 2.5 VS Probe - Hs-TERT
Probe Description	Target probe for Homo sapiens telomerase reverse transcriptase (TERT) transcript variant 1 mRNA. For Research Use Only.†	
Specificity	Homo sapiens TERT transcript variant 1 mRNA	
Species	Homo sapiens	
Target Gene	TERT	
Entrez Gene ID	7015	
Accession#	NM_198253.2	
Probe Region	164 - 3231	
ZZ Pair No.	20	

V-EBV-EBER1		
Catalog No.	310271	RNAscope Probe -V-EBV-EBER1
	310278	RNAscope 2.5 LS Probe - V-EBV-EBER1
	310279	RNAscope 2.5 VS Probe -V-EBV-EBER1
Probe Description	Target probe for Human herpesvirus 4 isolate SDTW400 (Epstein-Barr virus) encoded RNA-1 (EBER-1). For Research Use Only.† EBV target probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	EBV EBER-1 mRNA transcript	
Species	Epstein-Barr virus	
Target Gene	N/A	
Entrez Gene ID	N/A	
Accession#	KP195701.1	
Probe Region	39-120	
ZZ Pair No.	2	

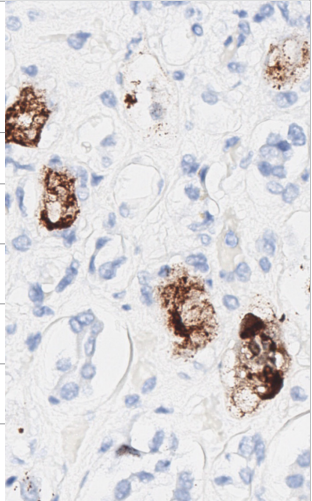
V-CMV		
Catalog No.	592701	RNAscope Probe - V-CMV
	592708	RNAscope 2.5 LS Probe - V-CMV
	592709	RNAscope 2.5 VS Probe - V-CMV
Probe Description	Target probe for Human cytomegalovirus Beta2.7 non-coding RNA transcript. For Research Use Only.† CMV probe is also available as an Analyte Specific Reagent (ASR) via Leica Biosystems.	
Specificity	CMV Beta2.7 non-coding RNA transcript	
Species	Human cytomegalovirus	
Target Gene	RNA2.7	
Entrez Gene ID	7755301	
Accession#	NC_006273.2	
Probe Region	4038-4158	
ZZ Pair No.	3	

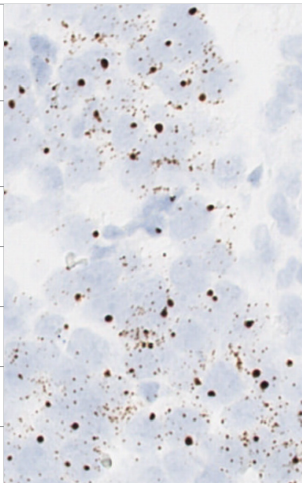
B-B.henselae - 23SrRNA		
Catalog No.	485621	RNAscope Probe - B-B. henselae-23SrRNA
	485628	RNAscope 2.5 LS Probe - B-B.henselae-23SrRNA
	485629	RNAscope 2.5 VS Probe - B-B.henselae-23SrRNA
Probe Description	Target probe for <i>Bartonella henselae</i> strain Houston-1 23S ribosomal RNA. For Research Use Only.†	
Specificity	<i>B. henselae</i> 23S ribosomal RNA	
Species	<i>Bartonella henselae</i> str. Houston-1	
Target Gene	23S rRNA	
Entrez Gene ID	N/A	
Accession#	NR_076339.1	
Probe Region	1210-1277	
ZZ Pair No.	1	

B-H.pylori - 16S		
Catalog No.	542931	RNAscope Probe - B-H. pylori-16S
	542938	RNAscope 2.5 LS Probe - B-H.pylori-16S
	542939	RNAscope 2.5 VS Probe - B-H.pylori-16S
Probe Description	Target probe for <i>Helicobacter pylori</i> strain ATCC 43504 16S ribosomal RNA. For Research Use Only.†	
Specificity	<i>Helicobacter pylori</i> ATCC 43504 16S ribosomal RNA	
Species	<i>Helicobacter pylori</i>	
Target Gene	16S rRNA	
Entrez Gene ID	N/A	
Accession#	NR_044761.1	
Probe Region	43-823	
ZZ Pair No.	2	

B-T.pallidum - 23SrRNA		
Catalog No.	579281	RNAscope Probe - B-T. pallidum-23SrRNA
	579288	RNAscope 2.5 LS Probe - B-T.pallidum-23SrRNA
	579289	RNAscope 2.5 VS Probe - B-T.pallidum-23SrRNA
Probe Description	Target probe for <i>Treponema pallidum</i> subsp. pallidum strain SS14 23S ribosomal RNA gene. For Research Use Only.†	
Specificity	<i>T. pallidum</i> subsp. pallidum strain SS14 23S rRNA	
Species	<i>T. pallidum</i> subsp. pallidum strain SS14	
Target Gene	23S rRNA	
Entrez Gene ID	NA	
Accession#	NR_076531.1	
Probe Region	121-228	
ZZ Pair No.	2	

V-RSV-NP		
Catalog No.	439861	RNAscope Probe - V-RSV-NP
	439868	RNAscope 2.5 LS Probe - V-RSV-NP
	439869	RNAscope 2.5 VS Probe - V-RSV-NP
Probe Description	Target probe for Respiratory syncytial virus isolate RSV Memphis-37 nucleoprotein (NP) mRNA. For Research Use Only.†	
Specificity	RSV NP mRNA	
Species	Respiratory Syncytial Virus	
Target Gene	NP	
Entrez Gene ID	N/A	
Accession#	KM360090.1	
Probe Region	8-1111	
ZZ Pair No.	20	

V-BKPyV-largeT		
Catalog No.	503521	RNAscope Probe - V-BK-PyV-largeT
	503528	RNAscope 2.5 LS Probe - V-BKPyV-largeT
	503529	RNAscope 2.5 VS Probe - V-BKPyV-largeT
Probe Description	Target probe for BK polyomavirus largeT mRNA transcript. For Research Use Only.†	
Specificity	BK polyomavirus largeT transcript	
Species	BK polyomavirus	
Target Gene	LargeT	
Entrez Gene ID	N/A	
Accession#	NC_001538.1	
Probe Region	48 - 2027	
ZZ Pair No.	28	

V-MCPyV-LT-ST-Ag-O1		
Catalog No.	833351	RNAscope Probe - V-MCPyV-LT-ST-Ag-O1
	833358	RNAscope 2.5 LS Probe - V-MCPyV-LT-ST-Ag-O1
	833359	RNAscope 2.5 VS Probe - V-MCPyV-LT-ST-Ag-O1
Probe Description	Target probe for Merkel cell polyomavirus isolate R17b. For Research Use Only.†	
Specificity	MCPyV isolate R17b LargeAg transcript	
Species	Merkel cell polyomavirus	
Target Gene	LargeAg	
Entrez Gene ID	N/A	
Accession#	NC_010277.2	
Probe Region	2-2445	
ZZ Pair No.	37	

bt

Still Need Antibodies? No Problem!

Primary Antibodies,
Secondary Antibodies,
and Isotype Controls
From R&D Systems



**View Our
Antibodies Here!**



Where Science Intersects Innovation™

Bio-Techne Brands

R&D Systems™ | Novus Biologicals™ | Tocris Bioscience™ | ProteinSimple™ | ACD™
bio-techne.com



Contact Us

Global info@bio-techne.com bio-techne.com/find-us/distributors
TEL +1 612 379 2956 North America TEL 800 343 7475
Europe | Middle East | Africa TEL +44 (0)1235 529449
China info.cn@bio-techne.com TEL +86 (21) 52380373

† For Research Use Only. Not for use in diagnostic procedures.

Trademarks and registered trademarks are the property of their respective owners.

MK 51-171 BR_ACD_AnatomicPathologyProbes_STRY0220208

bio-techne®