

RNAscope™ HiPlex Pro for COMET™ Ordering Checklist

This checklist provides a list of materials needed to execute the Multiomics protocol on COMET. The checklist notes materials required when RNAscope only as well as those needed to run RNA and protein.

Note: This checklist does not include materials needed for the preparation of the tissue blocks (sample fixation, embedding, sectioning). Tissue should be sectioned on Superfrost Plus slides.

Reagents/Materials Required from Advanced Cell Diagnostics

RNAscope HiPlex Pro for COMET 12-plex, 10-slide Reagent Kit (Cat. No 323680)

- | | |
|--|--------------------------------------|
| ☐ RNAscope HiPlex Pro for COMET 12-plex, 10-slide Detection Reagents | Cat. No. 323675 |
| ☐ RNAscope HiPlex Cleaving Stock Solution v2 (one ampoule required/run) | Cat. No. 324399 (1 box x 5 ampoules) |
| ☐ RNAscope HiPlex Pro Imaging Buffer (sufficient for 20-slides) | Cat. No 322080 |

Refer to page 2 for quantities provided in Reagent kit and storage information

Control Probes (Ready to Use, 1X), 10 slides

- | | |
|---|-----------------|
| ☐ RNAscope HiPlex12 CS Negative Control Probe | Cat. No. 324347 |
| ☐ RNAscope HiPlex12 CS Positive Control Probe – Hs | Cat. No. 324317 |
| ☐ RNAscope HiPlex12 CS Positive Control Probe – Mm | Cat. No. 324437 |
| ☐ RNAscope HiPlex12 CS Positive Control Probe – Rn | Cat. No. 324337 |

Target Probes (50X), 10 slides

- | | |
|---|------------------|
| ☐ RNAscope HiPlex CS Probe-[sp]-[target]-(T1...T12)
Example: RNAscope HiPlex CS Probe - Mm-Ifng-T10 | Cat. No. various |
|---|------------------|

Control Sample

- | | |
|---|--|
| ☐ RNAscope Control Slide -Human Hela Cell Pellet | Cat. No. 310045 (4, single-section slides) |
|---|--|

Requirements from Lunaphore

Instrument/Software (Required)

- | | | |
|--------------------------|---|--------------------------|
| ☐ | COMET™ 1.0 Instrument and control station | CM10-S |
| ☐ | COMET Software* | Version 1.6.0.0 or later |
| ☐ | PT Module (Epredia™) | AP07 |

*Note: license key necessary to unlock protocol, Contact Lunaphore

Consumables

- | | | |
|--------------------------|--|--------|
| ☐ | COMET chip | MK03 |
| ☐ | Multistaining Buffer (required) | BU06 |
| ☐ | Dewax and HIER Buffer H, pH 9 (Epredia™) | AR03 |
| ☐ | Imaging Buffer (required) | BU09 |
| ☐ | Elution Buffer Kit (required) | BU07-L |
| ☐ | Blocking Buffer Kit (optional) | BU10 |

Other Requirements for Multiomics on COMET

Reagents, consumables, equipment (required)

- | | | |
|--------------------------|--|---|
| ☐ | RNase Inhibitor, Murine | New England Biolabs (M0314S, M0314L) |
| ☐ | 20X SSC | Major Lab Supplier |
| ☐ | Xylene (or xylene-alternative, e.g. HistoClear) | Major Lab Supplier |
| ☐ | 100% EtOH | Major Lab Supplier |
| ☐ | Drying oven, capable of holding temperature at 60 +/- 1°C | Major Lab Supplier |
| ☐ | 50 mL conical tubes | See COMET user manual for compatible products |
| ☐ | 2 mL microtubes | See COMET user manual for compatible products |
| ☐ | Slide holders, staining dishes fume hood, etc. (for dewax) | Major Lab Supplier |
| ☐ | Tubes, bottles, reservoirs for buffer preparation | Major Lab Supplier |

Consumables (required when including protein cycles)

- | | | |
|--------------------------|----------------------|---------------|
| ☐ | Primary antibodies | User selected |
| ☐ | Secondary antibodies | User selected |

RNAscope HiPlex Pro for COMET Reagent Kit Components and Storage

Kit	Kit components	Quantity	Storage Temp.
RNAscope™ HiPlex Pro for COMET™ 12-plex, 10-slide Reagent Kit Cat. No. 323680	RNAscope™ HiPlex Pro for COMET™	1	2-8°C
	12-plex, 10-slide Detection Reagents Cat. No. 323675		
	RNAscope HiPlex Cleaving Stock Solution v2 Cat. No. 324399	2	15-30°C
	RNAscope HiPlex Pro Imaging Buffer Cat. No. 322080	1	15-30°C

Individual products with reagent information

Individual kits	Reagents	Volume/Amt.	Storage Temp.
RNAscope™ HiPlex Pro for COMET™	RNAscope™ HiPlex Pro AMP 1	5.2 mL x 1 bottle	2-8°C
	RNAscope™ HiPlex Pro AMP 2	5.2 mL x 1 bottle	2-8°C
	RNAscope™ HiPlex Pro AMP 3	5.2 mL x 1 bottle	2-8°C
12-plex, 10-slide Detection Reagents Cat. No. 323675	RNAscope™ HiPlex Pro Fluoro T1-T4	5.2 mL x 1 bottle	2-8°C
	RNAscope™ HiPlex Pro Fluoro T5-T8	5.2 mL x 1 bottle	2-8°C
	RNAscope™ HiPlex Pro Fluoro T9-T12	5.2 mL x 1 bottle	2-8°C
	RNAscope™ HiPlex PretreatPro™	5.2 mL x 1 bottle	2-8°C
	RNAscope™ HiPlex Pro Probe Diluent	5.2 mL x 1 bottle	2-8°C
	RNAscope™ post hybridization wash	8 mL x 1 bottle	2-8°C
	RNAscope™ HiPlex Pro DAPI (1 mg/mL)	1 mL x 1 bottle	2-8°C
RNAscope HiPlex Cleaving Stock Solution v2 Cat. No. 324399	5 Ampoules of cleaving solution	1 box	15-30°C
RNAscope HiPlex Pro Imaging Buffer Cat. No. 322080	RNAscope HiPlex Pro Imaging Buffer Part A	400 mg	15-30°C
	RNAscope HiPlex Pro Imaging Buffer Part B	100 mL	15-30°C