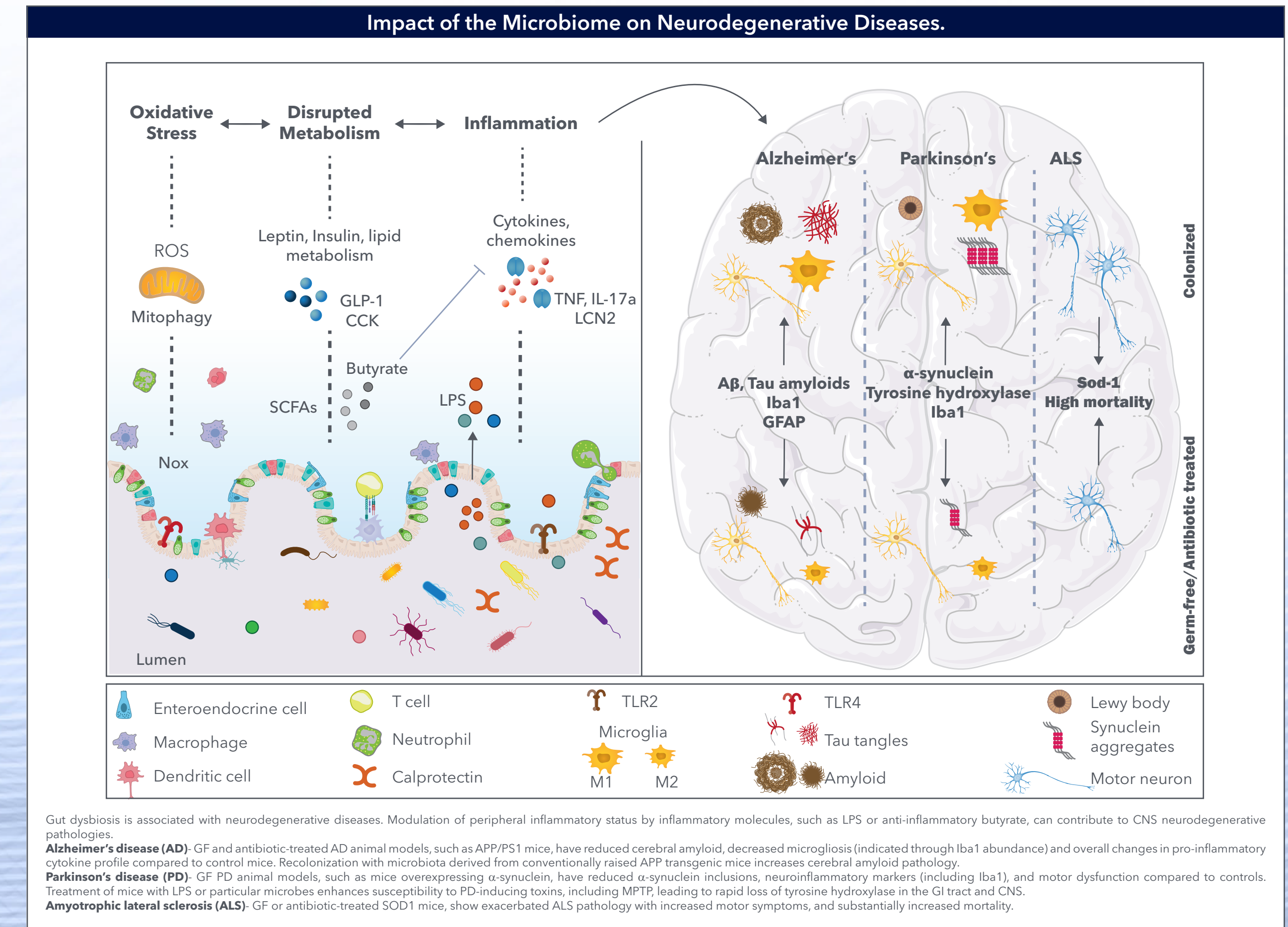
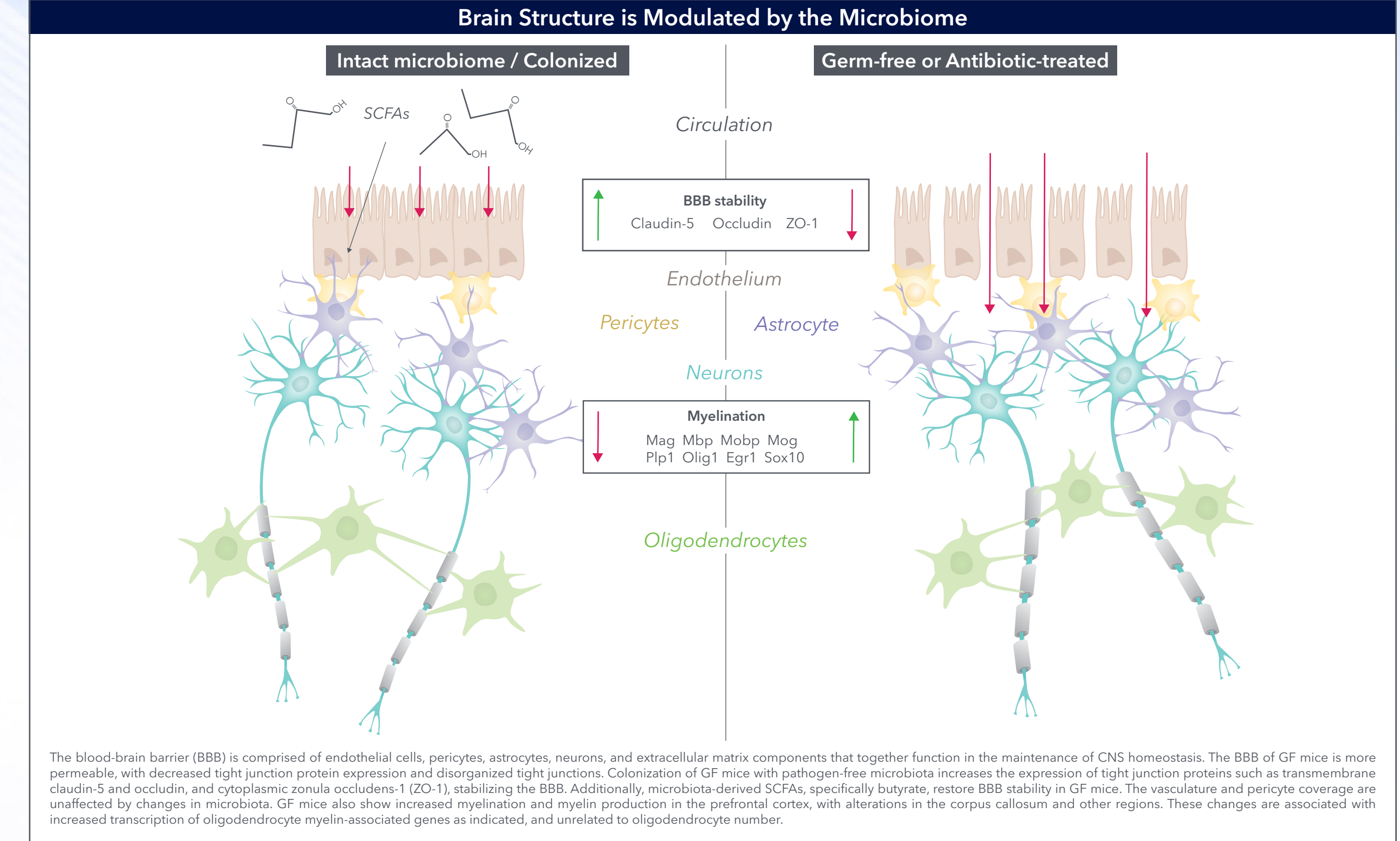
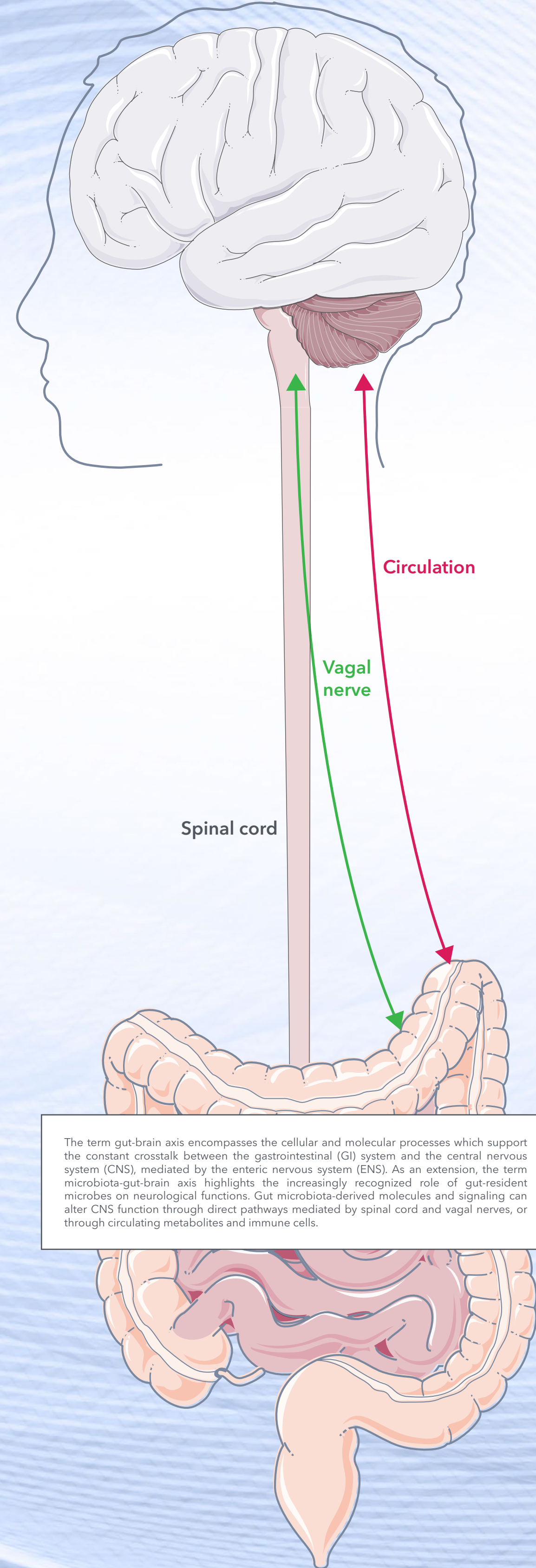
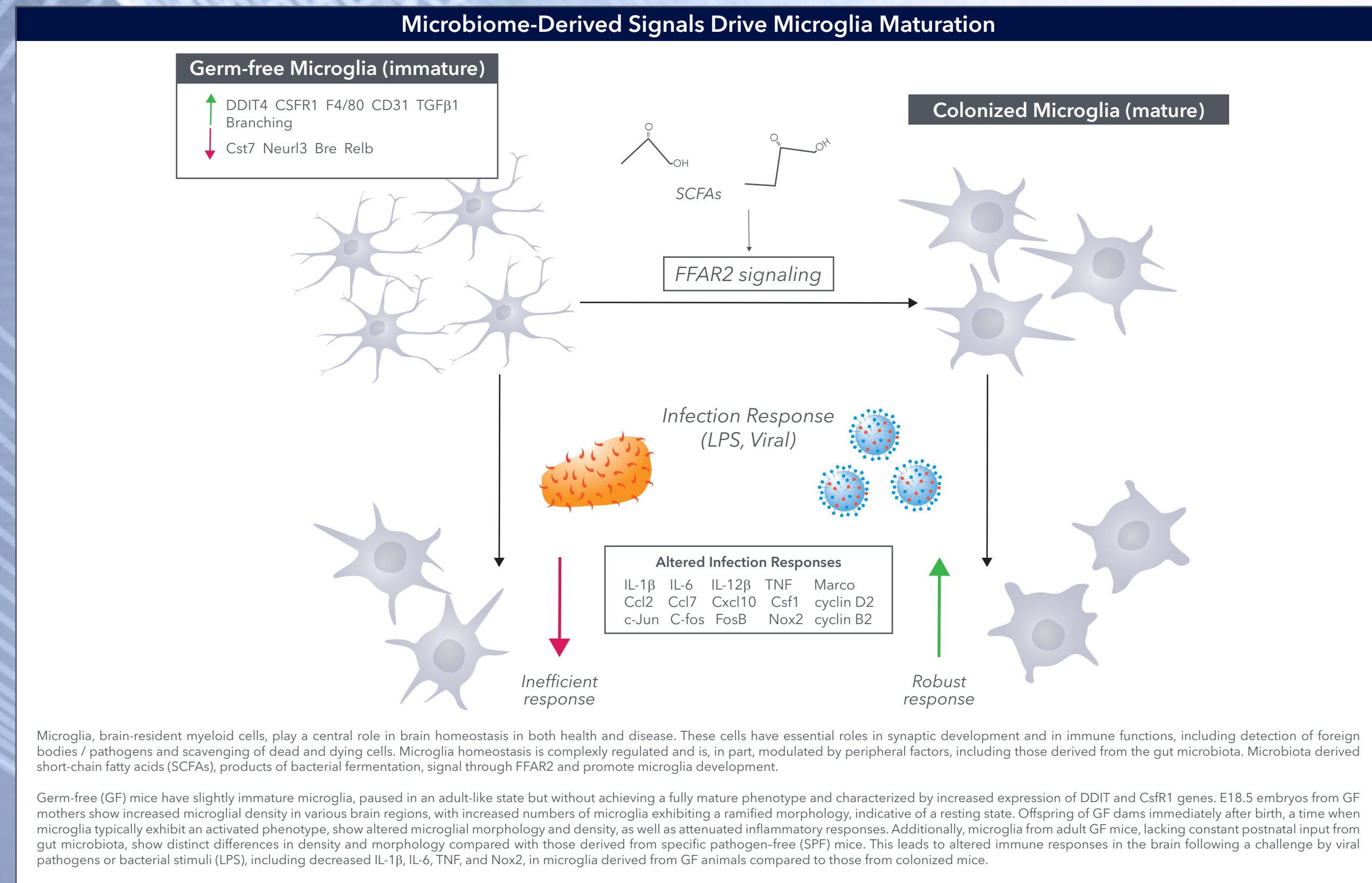
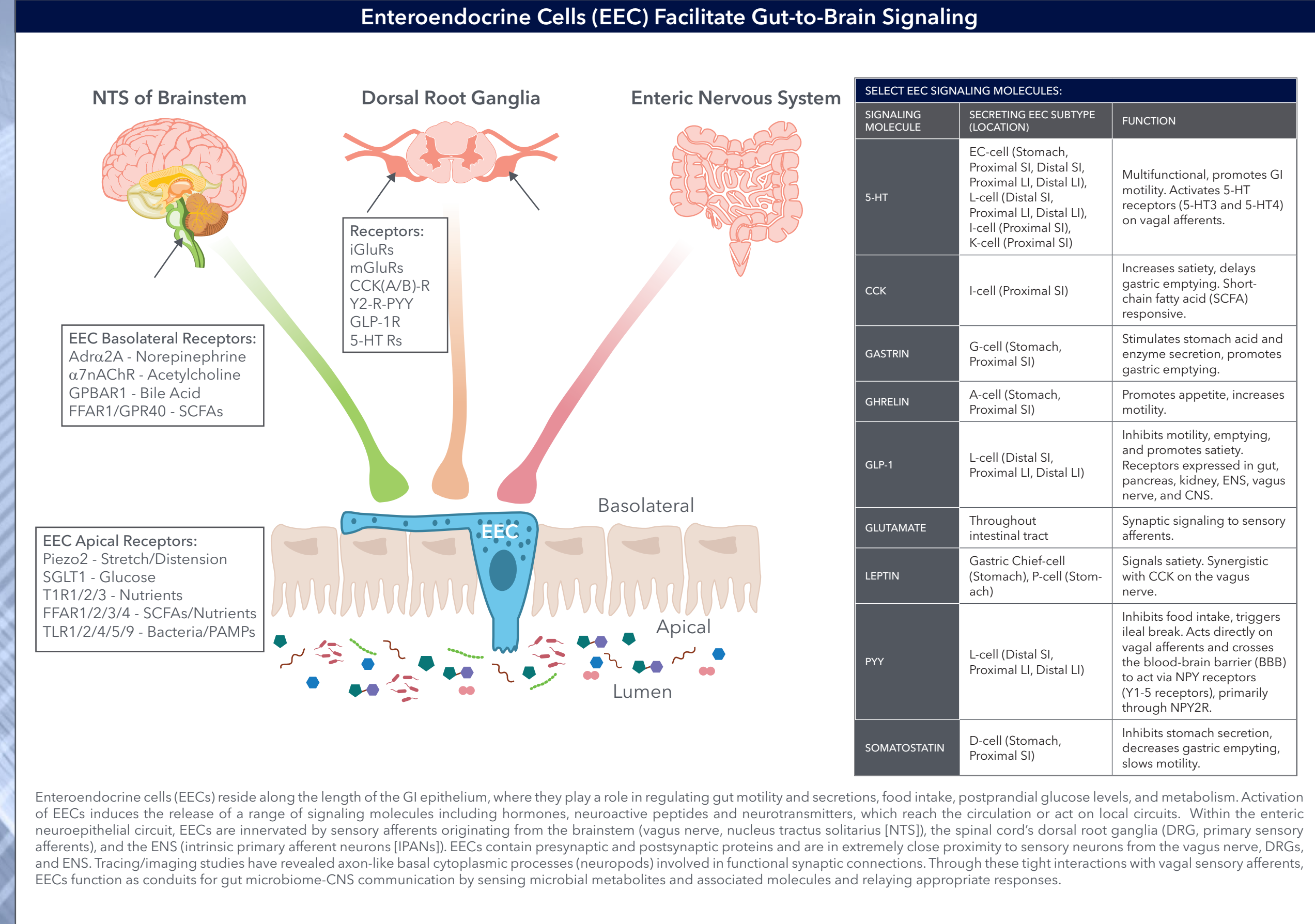


Timothy Sampson¹, Adam Hamilton², and Maria Elizabeth de Sousa Rodrigues³



NOTE: This poster conveys a general overview and should be considered neither comprehensive nor definitive. The details of this information are understood to be subject to interpretation.

ENTEROENDOCRINE CELLS (EEC) FACILITATE GUT-TO-BRAIN SIGNALING	
5-HT(R)	5-hydroxytryptamine (receptor)
α 7nAChR	Alpha-7 nicotinic receptor
Ad α 2A	Alpha-2A-adrenergic receptor
CCK(A/B)-R	Cholecystokinin (A/B) receptor
DRG	Dorsal root ganglia
EEC	Enteroendocrine cell
ENS	Enteric nervous system
FFAR1/2/3/4	Free fatty acid receptor 1/2/3/4
GI	Gastrointestinal
GLP-1(R)	Glucagon-like peptide 1 (receptor)
GPBAR1	G protein-coupled bile acid receptor 1
GPR40	G-protein-coupled receptor 40
iGluRs	Ionotropic glutamate receptors
IPAN	Intrinsic primary afferent neuron
LI	Large intestine
mGluRs	Metabotropic glutamate receptors
NPY (2R)	Neuropeptide Y (receptor Y2)
NTS	Nucleus tractus solitarius
PAMPs	Pathogen-associated molecular patterns
PYY	Peptide YY
SCFA	Short chain fatty acid
SGLT1	Sodium-glucose cotransporter 1
SI	Small intestine
T1R1/2/3	Taste receptor type 1, member 1/2/3
TLR1/2/4/5/9	Toll-like receptor 1/2/4/5/9
Y2-R - PYY	Y receptor type 2 - Peptide YY

MICROBIOME-DERIVED SIGNALS DRIVE MICROGLIA MATURATION	
Bre	Brain and reproductive organ-expressed
Ccl2	Chemokine (C-C motif) ligand 2
Ccl7	Chemokine (C-C motif) ligand 7
Csf1	Colony stimulating factor 1
CsfR1	Colony stimulating factor 1 receptor
CD31	Platelet endothelial cell adhesion molecule-1 (PECAM-1)
Cst7	Cystatin F
Cxcl10	C-X-C motif chemokine ligand 10
DDIT	DNA-damage inducible transcript
DDIT4	DNA-damage inducible transcript 4
F4/80	EGF-like module-containing mucin-like hormone receptor-like 1 (EMR1)
FFAR2	Free fatty acid receptor 2
FosB	FBJ murine osteosarcoma viral oncogene homolog B
GF	Germ-free
IL-1 β	Interleukin 1 beta
IL-6	Interleukin 6
IL-12 β	Interleukin 12 beta
LPS	Lipopolysaccharide
Marco	Macrophage receptor with collagenous structure
Neurl3	Neuralized E3 ubiquitin protein ligase 3
Nox2	NADPH oxidase 2
Relb	V-rel reticuloendotheliosis viral oncogene homolog B
SPF	Specific pathogen-free
TGF β 1	Transforming growth factor beta 1
TNF	Tumor necrosis factor

BRAIN STRUCTURE IS MODIFIED BY THE MICROBIOME	
BBB	Blood-brain barrier
CNS	Central nervous system
Egr1	Early growth response 1
GF	Germ-free
Mag	Myelin-associated glycoprotein
Mbp	Myelin basic protein
Mobp	Myelin-associated oligodendrocyte basic protein
Mog	Myelin oligodendrocyte glycoprotein
Olig1	Oligodendrocyte transcription factor 1
Plp1	Proteolipid protein 1
SCFA	Short-chain fatty acid
Sox10	SRY-related HMG-box 10
ZO-1	Zonula occludens-1

IMPACT OF THE MICROBIOME ON NEURODEGENERATIVE DISEASES	
A β	Beta-amyloid peptide
AD	Alzheimer's disease
ALS	Amyotrophic lateral sclerosis
APP	Amyloid precursor protein
CCK	Cholecystokinin
CNS	Central nervous system
GF	Germ-free
GFAP	Glial fibrillary acidic protein
GI	Gastrointestinal
GLP-1	Glucagon-like peptide 1
Iba1	Ionized calcium binding adapter molecule 1
IL-17a	Interleukin 17A
LCN2	Lipocalin-2
LPS	Lipopolysaccharide
MPTP	1-methyl-4-phenyl-1,2,3,6-tetrahydropyridine
Nox	NADPH oxidase
PD	Parkinson's disease
PS1	Presenilin 1
ROS	Reactive oxygen species
SCFA	Short-chain fatty acid
SOD1	Superoxide dismutase 1
TLR2	Toll-like receptor 2
TLR4	Toll-like receptor 4
TNF	Tumor necrosis factor