

# Cell-based assays for neurosciences



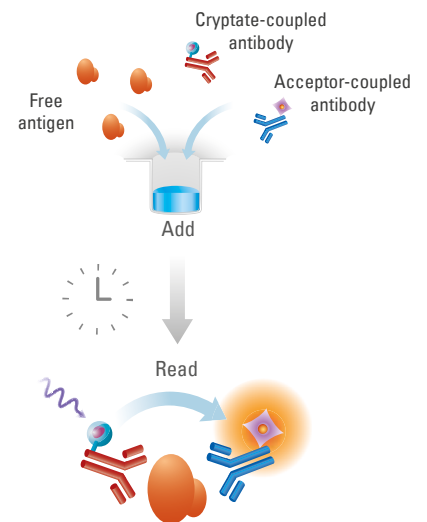
## HTRF® SOLUTIONS FOR NEUROSCIENCES

Whether you focus on Alzheimer's, Parkinson's, or Huntington's diseases, Friedreich's ataxia, other central nervous system (CNS) diseases, or disorders requiring pain management, Cisbio is committed to helping boost your research efforts.

Join the HTRF researcher community and move to HTRF®, a robust and reliable technology with a proven 20-year track record.

### Four reasons to choose Cisbio HTRF® solutions for your neuroscience projects:

- Over 110 cell-based neuroscience assays
- 10,000 HTRF® users worldwide and nearly 2,000 publications using HTRF®
- Dedicated scientist consultants to advise you and provide support
- High quality offer respecting same Quality Control process as Cisbio IVD products



### HTRF® technology, an effective alternative to ELISA & Western Blot

- Simple wash-free add-and-read protocol saves you time
- Low consumption preserves precious samples

NEUROSCIENCES

## OUR CNS PRODUCTS

### NEURODEGENERATIVE MARKERS

Alpha synuclein aggregation	CREB phospho-S133 & total	LRRK2 phospho-S935 & Total	Tau phospho-S202/T205
Apolipoprotein E (human)	Frataxin (human)	SMN (human)	Tau Aggregation
Amyloid $\beta$ 1-40 (human)	GSK3 alpha phospho-S21 & Total	Map Tau	
BACE1	GSK3 beta phospho-S9 & Total	Progranulin (human)	

### NEUROINFLAMMATION

CCL2/MCP1 (human & mouse)	IL1 $\beta$ (human & mouse)	IL12 (human)	TREM2 (human)
CCL4/MIP-1 $\beta$ (human)	IL2 (human & mouse)	IL17 (human & mouse)	
CCL20 (human)	IL4 (human)	IL18 (human)	
CXCL1/GRO $\alpha$ (human)	IL5 (human)	IL23 (human)	
G-CSF (human)	IL6 (human & mouse)	IL31 (human)	
IFN $\gamma$ (human & mouse)	IL10 (human)	TNF $\alpha$ (human, mouse & cynomolgus)	

### CELL SIGNALING

cAMP-Gs Dynamic	ATF2 phospho-T71	H2AX phospho-S139	mTOR phospho-S2448
cAMP-Gs HiRange	BAD phospho-S112	HER2 phospho-Y1221/1222 & Total	NF $\kappa$ B phospho-S536 & Total
cAMP-Gi	Beta Catenin phospho-T41/S33/37 & Total	Histone H3 phospho-T3 & Total	P38 phospho-T180/Y182
IP-One ELISA kit	cJun phospho-S63	Histone H3 phospho-S10	p70S6K phospho-T389
IP-One-Gq	CREB phospho-S133	IR beta phospho-Y1150/1151 & Total	SMAD3 phospho-S423/425 & Total
4EBP1 phospho-T37/46	EGFR phospho-Y1068, phospho-pan & Total	IRF3 phospho-S386	STAT1 phospho-Y701 & Total
ACC phospho-S79	EIF4E phospho-S209	IRS2 total	STAT3 phospho-Y705 & Total
AKT phospho-S473 & Total	Advanced phospho-ERK	IKKB phospho-S177/181 & Total	STAT5 phospho-Y694
AKT phospho-T308 & Total	ERK phospho-T202/Y204 & Total	JNK phospho-T183/Y185	Tau phospho-S202/T205
AKT1 phospho-S473 & Total	FGFR2 phospho-Y653/654 & Total	LRRK2 phospho-S935 & Total	ULK1 phospho-S758
AKT2 phospho-S473 & Total	GSK3 alpha phospho-S21 & Total	MEK1 phospho-S298, phospho-S218/222 & Total	VEGFR Phospho-Y1175
AKT3 phospho-S473 & Total	GSK3 beta phospho-S9 & Total	MEK1/2 phospho-S218/222	
AMPK alpha phospho-T172 & Total			

### NEUROTRANSMITTER RECEPTORS

Adenosine A2A binding	Cholecystokinin CCK1 binding	H3K4Me2 Cellular Assay	Muscarinic M5 binding
Adrenoceptor Beta 1 binding	Cholecystokinin CCK2 binding	Histamine H1 binding	Orexin OX1 binding
Adrenoceptor Beta 2 binding	Class A Orphans GPR1 binding	Histamine H3 binding	Orexin OX2 binding
Angiotensin AT2 binding	Dopamine D1 binding	Melanocortin MC4 binding	Serotonin 5HT1A binding
Bombesin BB1 binding	Dopamine D2 binding	Melanocortin MC5 binding	Serotonin 5HT1B binding
Bombesin BB2 binding	Dopamine D3a binding	Muscarinic M1 binding	Serotonin 5HT4 binding
Bradykinin B1 binding	Endothelin B binding	Muscarinic M2 binding	Tachykinin NK1 binding
Bradykinin B2 binding	Ghrelin GHSR1A binding	Muscarinic M4 binding	Vasopressin V2 binding

Discover all the details about Cisbio technology, services, and products for neurosciences on [neurosciences.cisbio.com](http://neurosciences.cisbio.com)

### EXPERT SERVICES TO SATISFY YOUR EXACT NEEDS



In addition to providing ready-to-use kits with all the necessary reagents, Cisbio's team of experienced experts works hand-in-hand with its clients to create custom solutions. When you have specific requirements for cellular phospho-proteins or other signaling proteins, Cisbio is the ideal trusted partner to develop the assays you need.

Keep your therapeutic research moving forward with faster, more robust, physiologically relevant results. Call us or visit [neurosciences.cisbio.com](http://neurosciences.cisbio.com)

#### FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 21 5018 9880 Japan +81-(0)43-306-8712 Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors

