

Cell-based assays for neurosciences



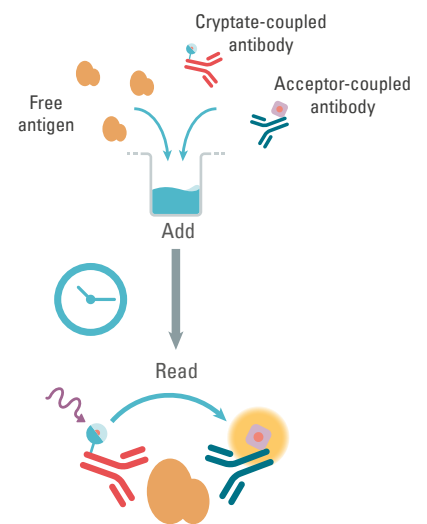
HTRF® SOLUTIONS FOR NEUROSCIENCES

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 | BACE1 | GSK3 beta phospho-S9 & Total | Progranulin (human) |
| Alpha Synuclein phospho-S129 & Total NEW | CREB phospho-S133 & total | LRRK2 phospho-S935 & Total | Tau phospho-S202/T205 NEW |
| Apolipoprotein E (human) | Frataxin (human) | SMN (human) | Tau phospho-S404 NEW |
| Amyloid β 1-40 (human) | GSK3 alpha phospho-S21 & Total | Map Tau | Tau Aggregation |

NEUROINFLAMMATION

| | | | |
|------------------------------|-----------------------------|--|---------------|
| CCL2/MCP1 (human & mouse) | IL1 β (human & mouse) | IL12 (human) | TREM2 (human) |
| CCL4/MIP-1 β (human) | IL2 (human & mouse) | IL17 (human & mouse) | |
| CCL20 (human) | IL4 (human) | IL18 (human) | |
| CXCL1/GRO α (human) | IL5 (human) | IL23 (human) | |
| G-CSF (human) | IL6 (human & mouse) | IL31 (human) | |
| IFN γ (human & mouse) | IL10 (human) | TNF α (human, mouse & cynomolgus) | |

CELL SIGNALING

| | | | |
|---|---|---|---|
| cAMP-Gs Dynamic | AMPK alpha phospho-T172 & Total | H2AX phospho-S139 | NFKB phospho-S536 & Total |
| cAMP-Gs HiRange | ATF2 phospho-T71 | HER2 phospho-Y1221/1222 & Total | P38 phospho-T180/Y182 |
| cAMP-Gi | BAD phospho-S112 | Histone H3 phospho-T3 & Total | p70S6K phospho-T389 |
| IP-One ELISA kit | Beta Catenin phospho-T41/S33/37 & Total | Histone H3 phospho-S10 | SMAD2 phospho-S465/467 & Total NEW |
| IP-One-Gq | cJun phospho-S63 | IR beta phospho-Y1150 /1151 & Total | SMAD3 phospho-S423/425 & Total |
| 4EBP1 phospho-T37/46 | CREB phospho-S133 | IRF3 phospho-S386 | STAT1 phospho-Y701 & Total |
| ACC phospho-S79 | EGFR phospho-Y1068, phospho-pan & Total | IRS2 total | STAT3 phospho-Y705 & Total |
| Alpha Synuclein phospho-S129 & Total NEW | EIF4E phospho-S209 | IKKB phospho-S177/181 & Total | STAT5 phospho-Y694 |
| AKT phospho-S473 & Total | Advanced phospho-ERK | JNK phospho-T183/Y185 | Tau phospho-S202/T205 |
| AKT phospho-T308 & Total | ERK phospho-T202/Y204 & Total | LRRK2 phospho-S935 & Total | Tau phospho-S404 NEW |
| AKT1 phospho-S473 & Total | FGFR2 phospho-Y653/654 & Total | MEK1 phospho-S298, phospho-S218/222 & Total | Tau phospho-S404 NEW |
| AKT2 phospho-S473 & Total | GSK3 alpha phospho-S21 & Total | MEK1/2 phospho-S218/222 | ULK1 phospho-S758 |
| AKT3 phospho-S473 & Total | GSK3 beta phospho-S9 & Total | mTOR phospho-S2448 | VEGFR Phospho-Y1175 |

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](https://www.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](https://www.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 to find a list of our regional distributors

