

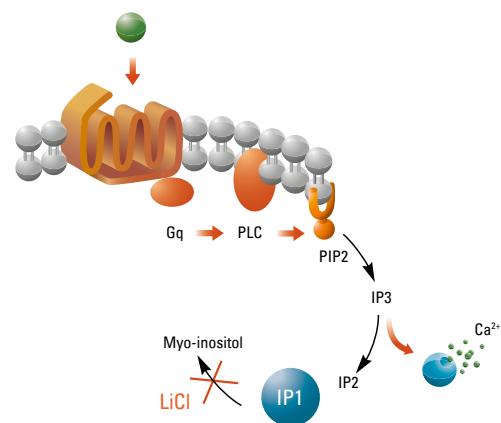
DON'T MISS A THING ON THE IP3 SIGNALING PATHWAY

HTRF® IP-One assays



IP-One is a unique assay to investigate Gq Protein-Coupled Receptors. IP-One assay detects inositol monophosphate, a stable downstream metabolite of IP3, induced by the activation of phospholipase C (PLC) cascade. Together with HTRF cAMP assays, IP-One provides a robust and reliable platform to assess GPCR targets, and a superior alternative to calcium sensing assays.

- Enables investigation of all classes of pharmaceutical compounds—agonists, antagonists, slow binders, allosteric modulators and inverse agonists
- Provides high hit confirmation—reduces false positives
- Can be used with a wide range of receptors and cell types, including endogenous receptors, transient or stable transfections
- Allows miniaturization to <math><10\ \mu\text{l}</math> for 1536 well protocols for HTS
- Also available in a traditional colorimetric ELISA format

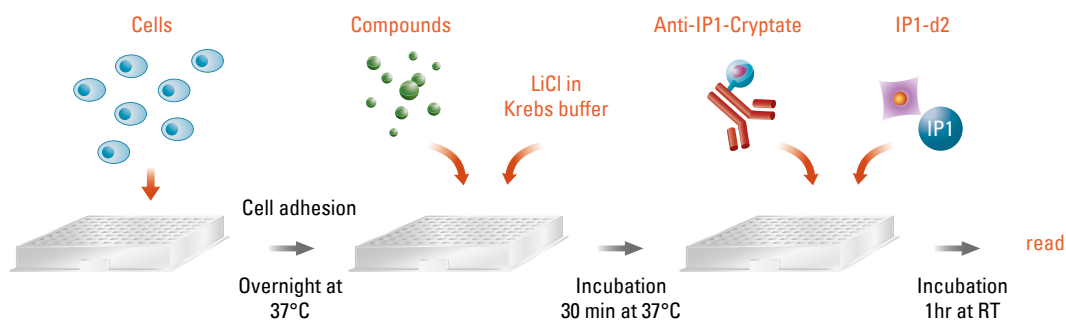


IP1 generation through the phospholipase C pathway.

A CELL-BASED, MIX AND READ ASSAY PROTOCOL

Upon stimulation of a Gq-coupled GPCR target, transduction signaling is triggered and the PLC pathway is activated. After lysis of the cell membrane, in the presence of LiCl preventing its degradation into myo-inositol, accumulation of inositol monophosphate (IP1) can be detected using the IP-One HTRF kit and most TR-FRET multimode plate readers.

Robust and efficient, the assay is based on a competitive format involving a specific antibody labeled with Terbium Cryptate (donor) and IP1 coupled to d2 (acceptor). The IP-One assay protocol consists of two incubation steps: cell stimulation by the ligand or target compounds, followed by IP1 detection using HTRF reagents. The assay can be run in a single microplate and requires only a single 1 hour incubation following cell stimulation.



ORDERING INFORMATION:

PRODUCT	TESTS	PART #	ADDITIONAL REAGENTS AVAILABLE	PART #
IP-One Tb kit	2 x 96	62IPAPEF	Lysis buffer 5 for IP-One Tb kit 1,000 tests—13 mL	62CL5FDD
IP-One Tb kit	1,000	62IPAPEB	Lysis buffer 6 for IP-One Tb kit 20,000 and 100,000 tests	62CL6FDF
IP-One Tb kit	20,000	62IPAPEC	IP-One calibrator	62IP1CDA
IP-One Tb kit	100,000	62IPAPEJ	IP-One stim. Buffer 5X—8 mL	62IP1FDC
IP-One Tb kit—10 packs 61PAPEJ	1,000,000	62IPAPEM	IP-One stim. Buffer 5X—100 mL	62IP1FDG
IP-One ELISA	96	72IP1PEA	IP-One Tb control	62IPATDA
IP-One ELISA	5 x 96	72IP1PED	96 half-well white plate cell culture—8 plates	66PL8ZZZ

For more information, please visit www.htrf.com/gpcr

INTERACTION IS EVERYTHING

From the chemistry of our assays to the collaborative nature of our relationship with customers, Cisbio is founded on the belief that “interaction is everything”. Talk to us. And discover a partner dedicated to providing exactly what you need:

- Convenient ready-to-use kits
- Custom assay capabilities
- Expert technical support
- Complete confidentiality
- Trusted and proven HTRF technology

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

