

## cAMP Hunter™ eXpress OPRL1 CHO-K1 GPCR Assay

**Catalog Number:** 95-0115E2

**Lot Number:** See Vial

**Contents:** 3.75 x 10<sup>6</sup> cells per vial in 0.2 mL

### Background

cAMP Hunter eXpress cells are engineered to overexpress naturally Gi or Gs coupled, wildtype GPCRs and are designed to detect changes in intracellular cAMP levels in response to agonist stimulation of the receptor. These cells have been modified to prevent long term propagation and expansion using a proprietary compound that has no apparent effect on assay performance.

### Product Information

**Target GPCR:** OPRL1

**Description:** Opioid receptor-like receptor

**Receptor Family:** Opioid

**Coupling:** Gi/Go

**Accession Number:** AY268428

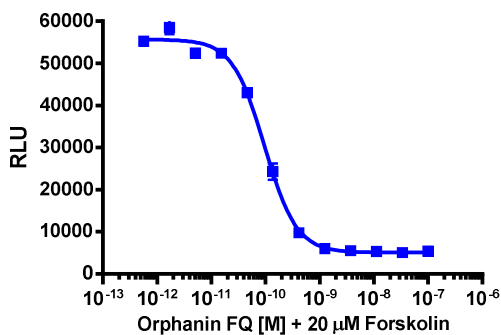
**GPCR Species:** Human

**Cell Type:** CHO-K1

**Storage:** Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

### Functional Performance

Cells were plated in a 96-well plate and stimulated with a control agonist (+ Forskolin for Gi targets), using the assay conditions described below. Following stimulation, signal was detected according to the recommended protocol. Please refer below for information on control compounds.



**Cell Number/Well:**

30000

**Control Agonist:**

Orphanin FQ

**Cell Plating Reagent:**

AssayComplete™ Cell Plating 2 Reagent

**Cell Incubation Time (Hours):**

24

**Agonist Incubation Time (minutes):**

30

**Agonist Incubation Temperature (°C):**

37

**EC<sub>50</sub> for Agonist Stimulation (nM):**

0.1

**Signal:Background at Agonist E<sub>max</sub>:**

10.2

### Additional Ligand Information

**Control Agonist:** Orphanin FQ

### Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit [discoverx.com/license](https://discoverx.com/license) for a list of products that are governed by limited use label license terms and relevant patent and trademark information.

**Ordering:** +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
**General product information:** [www.discoverx.com](https://www.discoverx.com)