

## cAMP Hunter™ eXpress AVPR2 CHO-K1 GPCR Assay

### MATERIALS PROVIDED

**Product Code:** 95-0150E2 **Lot Number:** See Vial

### BACKGROUND

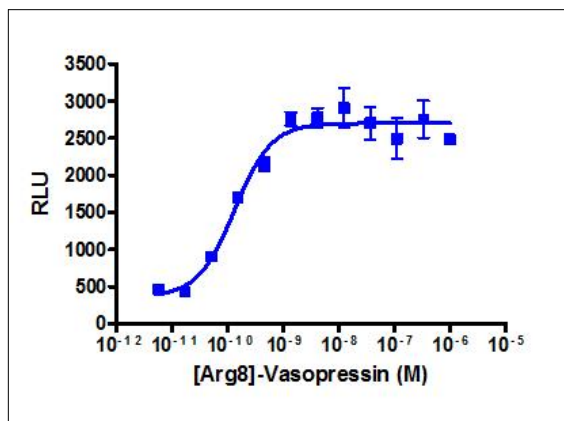
cAMP Hunter™ eXpress cells are engineered to overexpress naturally Gi or Gs coupled, wildtype GPCRs and are designed to detect changes in intracellular cAMP levels in response to agonist stimulation of the receptor. These cells have been modified to prevent long term propagation and expansion using a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

### PRODUCT INFORMATION

**Target GPCR:** AVPR2  
**Description:** Vasopressin receptor 2  
**Receptor Family:** Vasopressin and Oxytocin  
**Coupling:** Gs  
**Accession Number:** NM\_000054.4  
**GPCR Species:** Human  
**Cell Type:** CHO-K1  
**Format:** 1 vial, 2 vials or 10 vials, 3.75 x 10<sup>6</sup> cells per vial in 0.2 mL  
**Storage:** Short term (<2 weeks): Store at -80°C; Long term (>2 weeks): Store in vapor phase of liquid N<sub>2</sub>.

### 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:** [Arg<sup>8</sup>]-Vasopressin  
**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):** 0.13  
**Signal:Background at Agonist E<sub>max</sub>:** 5.4

### ADDITIONAL LIGAND INFORMATION

**Control Agonist:** [Arg<sup>8</sup>]-Vasopressin  
**Vendor:** DiscoverRx® (Part# 92-1105)

## LABEL LICENSE

For some products/cell lines, certain 3<sup>rd</sup> party gene specific patents may be required to use the cell line. It is the purchaser's responsibility to determine if such patents or other intellectual property rights are required.