

## MULTISCREEN™ STABLE CELL LINE RAT RECOMBINANT KAPPA OPIOID RECEPTOR

### PRODUCT INFORMATION

**Catalog Number:** C1352-1

**Lot Number:** C1352-1-061505

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO dhfr

**Transfection:** Full-length rat Oprk1 cDNA (GenBank Accession Number L22001)

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** Alpha-MEM, 10% FBS, 800 µg/mL G418

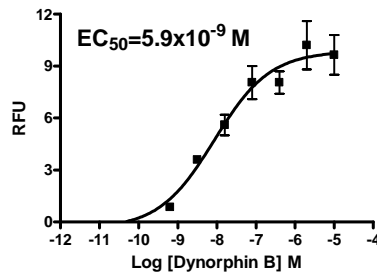
**Stability:** Stable after minimum of two months continuous growth

### Data sheet

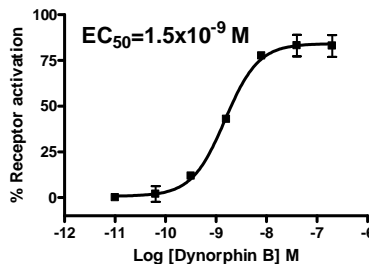
**Background:** Kappa opioid receptor (KOR) is a receptor for dynorphins. KOR inhibits neurotransmitter release by reducing calcium ion currents and increasing potassium ion conductance and may play a role in arousal and regulation of autonomic and neuroendocrine functions. Some studies suggest that the stimulation of kappa opioid receptors improves memory dysfunctions resulting from the blockade of muscarinic M1 receptors. In addition, there is evidence that kappa opioid receptor agonists attenuate several behavioral responses induced by drugs of abuse, raising the possibility that kappa opioid receptor agonists may be useful for the treatment of dependence on drugs of abuse.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC).

### References:

Ukai *et al.* (1995) Kappa-Opioid receptor agonists improve pirenzepine-induced disturbance of spontaneous alternation performance in the mouse. *Eur J Pharmacol* 281:173-178.

Hasebe *et al.* (2004) Possible pharmacotherapy of the opioid kappa receptor agonist for drug dependence. *Ann N Y Acad Sci* 1025:404-413.

**FOR RESEARCH USE ONLY.**

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com  
sales@multispaninc.com  
support@multispaninc.com

Ver. June 2005

Phone: +1 (510) 887-0817  
Fax: +1 (510) 887-0863  
26219 Eden Landing Road  
Hayward, CA 94545-3718  
U.S.A.