

**MULTISCREEN™ STABLE CELL LINE**  
**RAT RECOMBINANT B2 RECEPTOR**

**Data Sheet**

**PRODUCT INFORMATION**

**Catalog Number:** Cr1199

**Lot Number:** Cr1199-031516

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

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

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length rat B2 cDNA (GenBank Accession Number NM\_173100.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

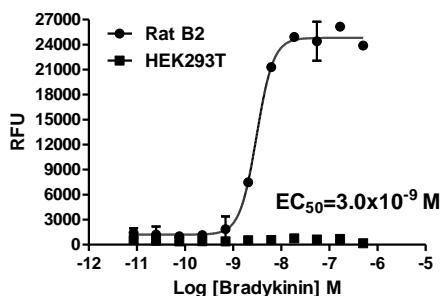
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

**Stability:** Stability in progress

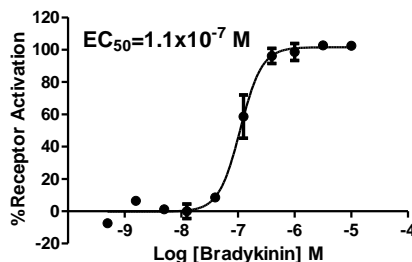
**Background:** Bradykinin receptor B2 is a constitutively expressed G protein-coupled receptor. In rats, the B2 receptors is present on non-peptidergic C-neurons, peptidergic C-neurons, dendrites and axons in A-neurons, NGF-responsive nociceptors. Recent research on rat TG neurons revealed that the B2 receptor is essential in the early stages of general pain generation and that endogenous bradykinin preferentially activates the B2 receptor over the B1 receptor. The activation of the B2 receptor by bradykinin stimulates neurogenic chloride secretion in the rat colon through COX increasing PGE2 production.

**Application:** Functional assays

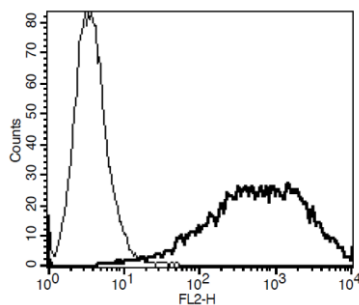
**Figure 1**



**Figure 2**



**Figure 3**



**Figure 1.** Dose-dependent calcium flux upon treatment with ligand, monitored with FLIPR.  
**Figure 2.** Dose-dependent increase of intracellular cAMP level upon treatment with ligand.  
**Figure 3.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

**References:**

Kawaguchi, A., Sato, M., Kimura, M., Yamazaki, T., Yamamoto, H., Tazaki, M., ... Shibukawa, Y. (2015). Functional expression of bradykinin B<sub>1</sub> and B<sub>2</sub> receptors in neonatal rat trigeminal ganglion neurons. *Frontiers in Cellular Neuroscience*, 9, 229. <http://doi.org/10.3389/fncel.2015.00229>

Radka Zubakova, Andreas Gille, Alexander Faussner, Ulrich Hilgenfeldt. (2008). Ca<sup>2+</sup> signalling of kinins in cells expressing rat, mouse and human B<sub>1</sub>/B<sub>2</sub>-receptor, *International Immunopharmacology*, 8(2), 276-281.

Qu, M.-H., Ji, W.-S., Zhao, T.-K., Fang, C.-Y., Mao, S.-M., & Gao, Z.-Q. (2016). Neurophysiological mechanisms of bradykinin-evoked mucosal chloride secretion in guinea pig small intestine. *World Journal of Gastrointestinal Pathophysiology*, 7(1), 150–159.

**FOR RESEARCH USE ONLY.**