

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT B1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1198a

**Lot Number:** C1198a-061510

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

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

**Host cell:** HEK293T

**Transfection:** Full-length Human BDKRB1 cDNA (GenBank Accession Number NM\_000710) with FLAG-tag sequence at the N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

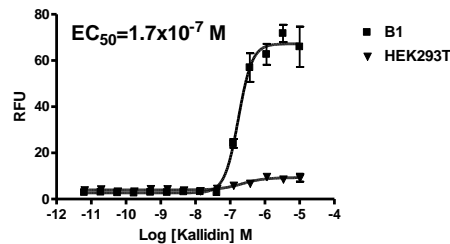
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

**Stability:** Stable in culture for minimum of two months

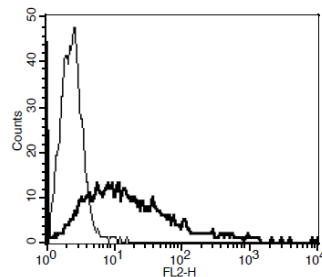
**Background:** Bradykinin B1 is a receptor for bradykinin and mediates responses to pathophysiological conditions such as inflammation, trauma, burns, shock, and allergy. Receptors for B1 are normally absent, but become highly upregulated after inflammatory stimuli. B1 receptor agonists may have important clinical value in the treatment of chronic pain and inflammatory disorders.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** 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:

- Bastian *et al.* (1997) Stable expression of human kinin B1 receptor in 293 cells: pharmacological and functional characterization. *Br J Pharmacol* 122:393-399.
- Calixto *et al.* (2004) Kinin B1 receptors: Key G-protein-coupled receptors and their role in inflammatory and painful processes. *Br J Pharmacol* 143:803-818.
- Menke *et al.* (1994) Expression cloning of a human B1 bradykinin receptor. *J Biol Chem* 269:21583-21586.

**FOR RESEARCH USE ONLY.**

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