

MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT B1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: Cm1198

Lot Number: Cm1198-082313

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Full-length Mouse BDKRB1 cDNA (GenBank Accession Number NM_007539.2) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin

Stability: In progress

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

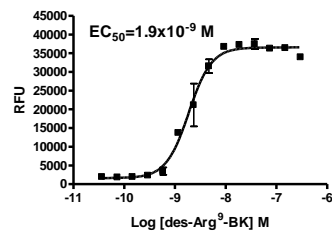


Figure 2

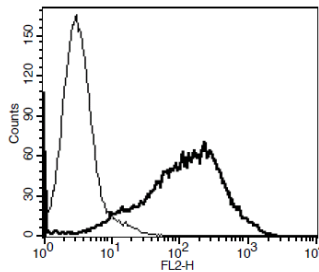


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

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.

Li *et al.* (1998) Involvement of bradykinin B₁ and B₂ receptors in relaxation of mouse isolated trachea. *Br J Pharmacol* 123:1337-1342.

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.