

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT B2 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DH1199

Lot Number: DH1199-100515

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Full-length Human BDKRB2 cDNA (GenBank Accession Number NM_000623) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable for 1-2 days after thawing

Background: Bradykinin receptor B2 is a G protein-coupled receptor for bradykinin. B2 receptor agonists may have important clinical value in the treatment and prevention of various cardiovascular disorders such as hypertension, ischemic heart disease, left ventricular hypertrophy, ventricular remodeling and congestive heart failure, as well as diabetic disorders by mimicking the reported beneficial effects of bradykinin. Blocking bradykinin B2 receptors after experimental cerebral ischemia reduces brain edema, infarct volume and neuronal necrosis, and improves neurological outcome. Thus, B2 antagonists may be a promising new class of compounds for clinical use after the onset of cerebral ischemia.

Application: Functional assays

Figure 1

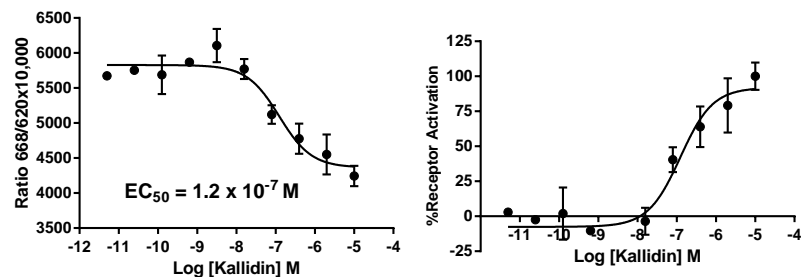


Figure 1. Dose-dependent stimulation of intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Hess *et al.* (1992) Cloning and pharmacological characterization of a human bradykinin (BK-2) receptor. *Biochem Biophys Res Commun* 184:260-268.

Heitsch (2003) The therapeutic potential of bradykinin B2 receptor agonists in the treatment of cardiovascular disease. *Expert Opin Investig Drugs* 12:759-770.

Sobey (2003) Bradykinin B2 receptor antagonism: a new direction for acute stroke therapy? *Br J Pharmacol* 139:1369-1371.

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.