

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT NK2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1304

Lot Number: C1304-073005

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human TACR2 cDNA (GenBank Accession Number NM_001057) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable after minimum of two months continuous growth

Background: The tachykinin receptor 2 (TACR2 or NK2) is a receptor for the tachykinin family of peptide neurotransmitters. NK2 receptor selectively binds neurokinin A (substance K), which is predominating in human and animal airways. It has been shown to regulate circular muscle contraction. NK2 can be found in tracheal and gastric tissues.

Application: Functional assays

Figure 1

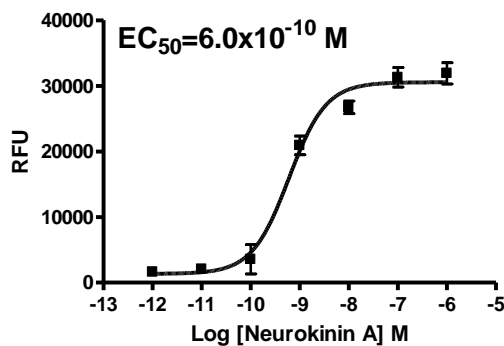


Figure 2

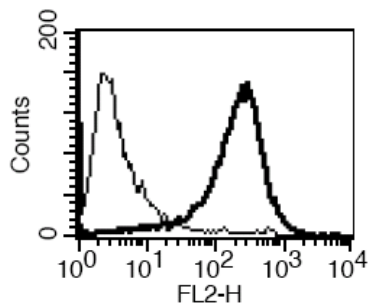


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick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

References:

Sasai and Nakanishi (1989) Molecular characterization of rat substance K receptor and its mRNAs. *Biochem Biophys Res Commun* 165:695-702.

Morten *et al.* (1991) Human NK-2 receptor gene maps to chromosome region 10q11-21. *Hum Genet* 88:200-203.

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.