

MULTISCREEN™ STABLE CELL LINE RAT RECOMBINANT NK3 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: Cr1305

Lot Number: Cr1305-011122

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: CellBanker2

Host cell: HEK293T

Transfection: Full-length Rat NK3 cDNA (GenBank Accession Number NM_017053.1) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: In progress

Background: Three membrane receptors that recognize these tachykinins have been described, termed TACR1, TACR2, and TACR3 (or NK3). Their endogenous ligands are substance P, neurokinin A, and neurokinin B, respectively. Interest in the tachykinins and their respective receptors has been stimulated by their potential role in bronchial asthma.

Application: Functional assays

Figure 1.

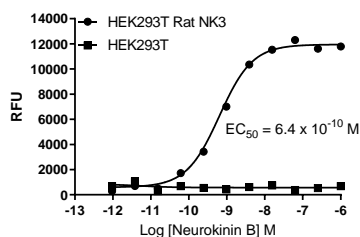


Figure 2.

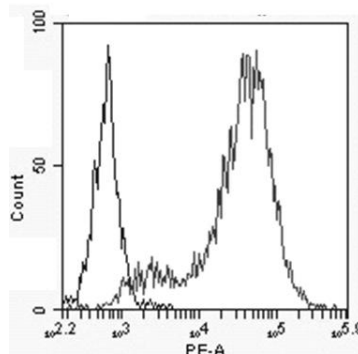


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Black line: parental cells; gray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

References:

Hopkins *et al.* (1991) Isolation and characterisation of the human lung NK-1 receptor cDNA. *Biochem Biophys Res Commun* 180:1110-1117.

Huang *et al.* (1992) cDNA sequence and heterologous expression of the human neurokinin-3 receptor. *Biochem Biophys Res Commun* 184:966-972.

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.