

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT NK3 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1305

Lot Number: MC1305-052025

Quantity: 1 vial (16.01 mg/mL; in 1mg, 5mg, or 10 mg)

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, 25 mM Sucrose (pH 7.2)

Host cell: HEK293T

Transfection: Expression vector containing full-length human TACR3 cDNA (GenBank Accession Number NM_001059.1) with FLAG tag sequence at N-terminus

Recommended Storage: -80°C upon receiving. Avoid repeated freeze-thaw.

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

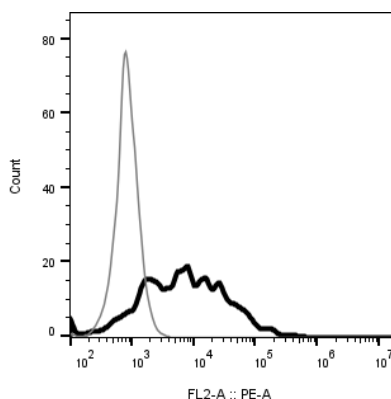


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

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.