

## MULTISCREEN™ DIVISION ARRESTED CELL LINE HUMAN RECOMBINANT NK2 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** DC1304

**Lot Number:** DC1304-020422

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cellbanker 2  
(Amsbio)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human TACR2 cDNA (GenBank Accession Number NM\_001057) with FLAG tag sequence at N-terminus

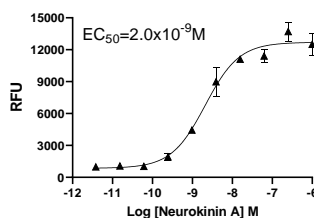
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS

**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 1.** Dose-dependent calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

#### 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.