

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT TRKA RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CK1001-1b

Lot Number: CK1001-1b-102615

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human FGFR3 cDNA (GenBank Accession Number NM_001012331) and vector pGL4.30[luc2P/NFAT-RE/Hygro] (GenBank Accession Number DQ904462)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 400 $\mu\text{g}/\text{mL}$ G418; 250 $\mu\text{g}/\text{mL}$ hygromycin

Stability: In progress

Background: Neurotrophic tyrosine kinase receptor type 1 (also called NTRK1; MTC; p140-TrkA; TRK; Trk-A; TRK1; TRKA) is a member of the neurotrophic tyrosine kinase receptor family. NTRK1 is a membrane bound receptor that phosphorylates itself and members of the MAPK pathway upon neurotrophin binding. Genetic variations in NTRK1 have been linked to anhidrosis, congenital insensitivity to pain, cancer, and mental retardation.

Application: Functional assay

Figure 1

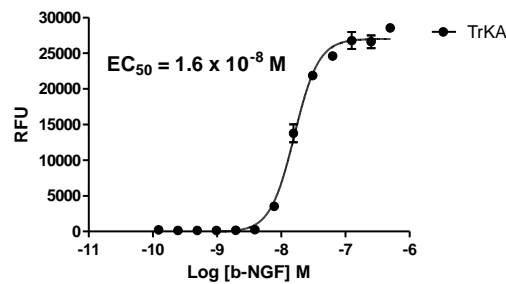


Figure 1. Dose-dependent calcium flux upon treatment with ligand, monitored with FLIPR.

References:

Bradshaw, R.A. et al. "Receptor Tyrosine Kinase Signaling Mechanisms: Devolving TrkA Responses with Phosphoproteomics." *Advances in biological regulation* 53.1 (2013): 87–96. *PMC*. Web. 4 Feb. 2015.

Mardy, S et al. "Congenital Insensitivity to Pain with Anhidrosis: Novel Mutations in the TRKA (NTRK1) Gene Encoding a High-Affinity Receptor for Nerve Growth Factor." *American Journal of Human Genetics* 64.6 (1999): 1570–1579. Print.

Marlin, M. Caleb, and Guangpu Li. "Biogenesis and Function of the NGF/TrkA Signaling Endosome." *International review of cell and molecular biology* 314 (2015): 239–257. *PMC*. Web. 4 Feb. 2015.

FOR RESEARCH USE ONLY.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.