

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT FGFR3 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CK1002-1

Lot Number: CK1002-1-020615

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_000142) 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: Fibroblast growth factor receptor 3 (FGFR3) is a member of the fibroblast growth receptor family. FGFR3 is a receptor tyrosine kinase that is activated by both acidic and basic fibroblast growth factors. FGFR3 plays an important role in bone maintenance and development. Mutations in FGFR3 have been linked to osteochondrodysplasia, craniosynostosis, and achondroplasia. Recent studies have also shown that somatic mutations in the FGFR3 gene are associated with bladder cancer, cervical cancer, and multiple myeloma.

Application: Functional assay

Figure 1

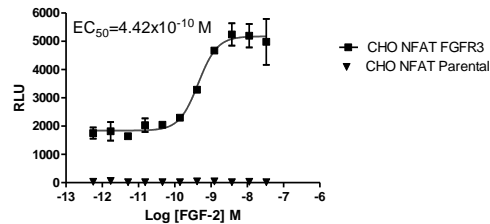
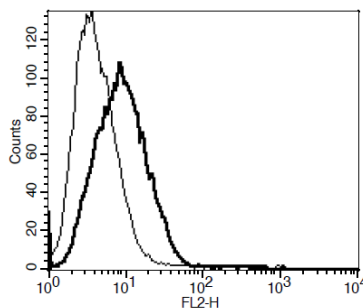


Figure 1. Ligand-induced dose-dependent luciferase activity measured by FlexStation III after 24 hours of treatment with the ligand in starvation media containing 0.2% FBS. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FGFR3 antibody. Thin line: parental cells; thick line: receptor-expressing cells.

Figure 2



References:

- Foldynova-Trantirkova, Silvie, William R. Wilcox, and Pavel Krejci. "Sixteen Years and Counting: The Current Understanding of Fibroblast Growth Factor Receptor 3 (FGFR3) Signaling in Skeletal Dysplasias." *Human mutation* 33.1 (2012): 29–41. *PMC*. Web. 3 Feb. 2015.
- Keegan, K et al. "Isolation of an Additional Member of the Fibroblast Growth Factor Receptor Family, FGFR-3." *Proceedings of the National Academy of Sciences of the United States of America* 88.4 (1991): 1095–1099. Print.
- Qing, Jing et al. "Antibody-Based Targeting of FGFR3 in Bladder Carcinoma and T(4;14)-Positive Multiple Myeloma in Mice." *The Journal of Clinical Investigation* 119.5 (2009): 1216–1229. *PMC*. Web. 3 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.