

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT TP RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1365

Lot Number: C1365-111909

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human TBXA2R cDNA (GenBank Accession Number NM_001060.4) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable in culture for minimum of two months

Background: The human TP (thromboxane A₂) receptor is a very potent stimulator of platelet aggregation and a constrictor of vascular and respiratory smooth muscles. It has been shown to be a mediator in diseases such as myocardial infarction, stroke and bronchial asthma. TP receptors can be found on platelets, as well as macrophages, monocytes, vascular endothelial cells, and smooth muscle cells. TP receptor antagonists may also play a role in the treatment of atherothrombosis and stroke prevention.

Application: Functional assays

Figure 1

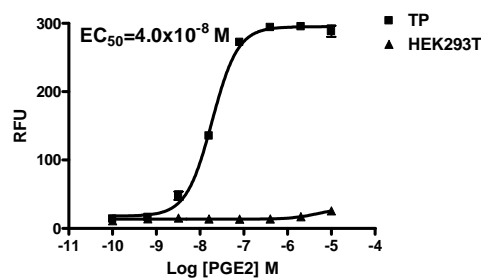


Figure 2

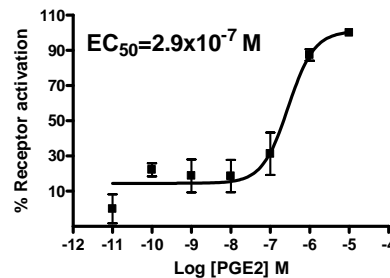


Figure 3

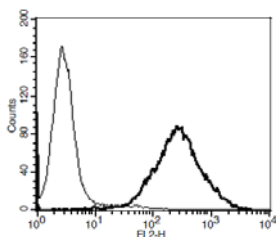


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent accumulation of intracellular IP1 upon treatment with ligand, measured with IP-one Tb kit (Cisbio 62IPAPC). **Figure 3.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Chamorro (2009) TP receptor antagonism: a new concept in atherothrombosis and stroke prevention. *Cerebrovasc Dis* 27 Suppl 3:20-27.

Hirata *et al.* (1991) Cloning and expression of cDNA for a human thromboxane receptor. *Nature* 349:617-20.

Nusing *et al.* (1993) Characterization and chromosomal mapping of the human thromboxane A₂ receptor gene. *J Biol Chem* 268:25253-25259.

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.