

**MULTISCREEN™ STABLE CELL LINE
RAT RECOMBINANT ET_A RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: Cr1216-1

Lot Number: Cr1216-1-083112

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Full-length Rat EDNRA cDNA (GenBank Accession Number NM_012550.2) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DME/F12, 10% FBS, 10 µg/mL puromycin

Stability: In progress

Background: Endothelin is a hormone produced predominantly by endothelial cells that have been recognized to play a significant role in the development of several cardiovascular disease states. Its receptor ET_A has been reported in a variety of tissues including placenta, uterus, testis, heart, adrenal gland, lung, kidney, ovary, skin, liver, bladder, large blood vessels, cerebellum, neurons, and meningioma tumor cells. In the vasculature, ET_A receptors residing on smooth muscle cells mediate prolonged vasoconstriction. Furthermore, activation of the ET_A mediates events that regulate mitogenesis, apoptosis, angiogenesis and metastasis in tumors.

Application: Functional assays

Figure 1

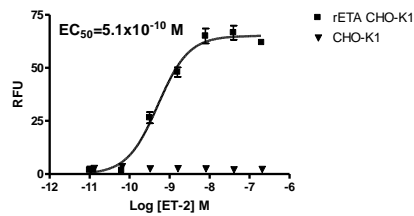


Figure 2

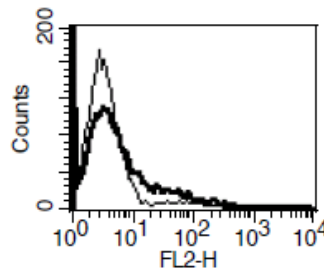


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Dupuis (2000) Endothelin receptor antagonists and their developing role in cardiovascular therapeutics. *Can J Cardiol* 16:903-910.

Wu-Wong (2002) Endothelin receptor antagonists as therapeutic agents for cancer. *Curr Opin Investig Drugs* 3:1234-1239.

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.