

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT SST5 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1349-1

**Lot Number:** DC1349-1-110218

**Quantity:** 1 vial ( $4 \times 10^6$ ) frozen cells

**Freeze Medium:** Cellbanker 2

**Host cell:** CHO-K1

**Transfection:** Expression vector containing full-length human SSTR5 cDNA (GenBank Accession Number NM\_001053) with FLAG tag sequence at N-terminus

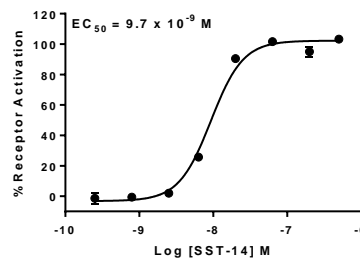
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM/F12, 10% FBS

**Background:** Somatostatin receptors (SSTRs) are activated by somatostatin secreted from nerve and endocrine cells. SSTRs are expressed in a tissue-specific manner and involved in the regulation of secretion of insulin, glucagon and growth hormone as well as cell growth induced by neuronal excitation in both the central and peripheral nervous systems. Aberrant expression of somatostatin receptors is known in a large number of human tumors. SSTR5 is expressed in adult pituitary gland, heart, small intestine, adrenal gland, cerebellum and fetal hypothalamus. SSTR5 is a candidate gene for bipolar affective disorder as well as for other neuropsychiatric disorders. Expression of SSTR5 may be important in the growth inhibitory effect of somatostatin in human pancreatic cancer.

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

**References:**

Ardjomand *et al.* (2003) Expression of somatostatin receptors in uveal melanomas. *Invest Ophthalmol Vis Sci* 44:980-987.

Ballare *et al.* (2001) Mutation of somatostatin receptor type 5 in an acromegalic patient resistant to somatostatin analog treatment. *J Clin Endocr Metab* 86:3809-3814.

**FOR RESEARCH USE ONLY.**

© 2018 by Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.