

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT SST3 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CG1347-1

Lot Number: CG1347-1-052715

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1 Gα16

Transfection: Expression vector containing full-length human SSTR3 cDNA (GenBank Accession Number: AY322541) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: Stable for 1-2 days after thawing

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. SSTR3 mRNA is detected in the brain and pancreatic islets. Aberrant expression of somatostatin receptors is known in a large number of human tumors.

Application: Functional assay

Figure 1

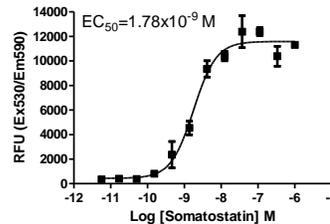


Figure 2

Figure 1. Dose-dependent stimulation of intercellular calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

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

Brinkmeier and Camper (1997) Localization of somatostatin receptor genes on mouse chromosomes 2, 11, 12, 15, and 17: correlation with growth QTLs. *Genomics* 43: 9-14.

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.