

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT CCK1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1038-1

Lot Number: C1038-1-120810

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Full-length human CCKAR cDNA (GenBank accession Number NM_000730.1) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 10 μ g/mL puromycin

Stability: In progress

Background: Cholecystikinin (CCK) is a regulatory peptide hormone with many functions in the gastrointestinal tract and brain. CCK1 (also known as CCKA) receptors are present in the gall bladder, gastric smooth muscle cells and mucosa, myenteric neurons, and in the brain. CCK regulates gastric acid secretion, gastric emptying and gall bladder motility in the gastrointestinal tract. Clinical studies show that CCK1 may be involved in many types of diseases, such as pancreatic disorders, motility disorders, tumor growth and regulation of satiety.

Application: Functional assays

Figure 1

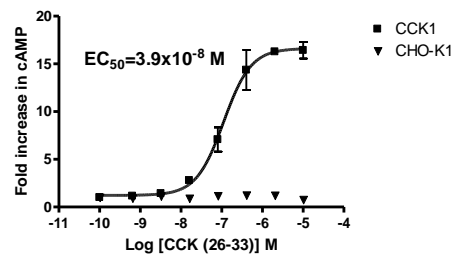


Figure 2

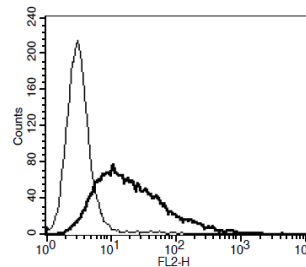


Figure 1. Dose-dependent inhibition of intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 2.** Receptor Expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Berna, MJ *et al.* (2007) Role of CCK/gastrin receptors in gastrointestinal/metabolic diseases and results of human studies using gastrin/CCK receptor agonists/antagonists in these diseases. *Curr Top Med Chem* 7(12):1211-31.

de Weerth, A *et al.* (1993) Molecular cloning, functional expression and chromosomal localization of the human cholecystikinin type A receptor. *Biochem Biophys Res Commun* 194(2):811-8.

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.