

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT OX2 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1034-1

Lot Number: 03/12/13

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Full-length Human HCRTR2 cDNA (GenBank Accession Number NM_001526.2) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DME/F12, 10% FBS

Stability: Stable in culture for 1-2 days after thawing

Background: Orexins, or hypocretins, produced by a small group of neurons in lateral hypothalamus, are involved in many physiological functions such as feeding, energy metabolism, sleep, arousal, reward, substance abuse, stress, as well as sympathetic and cardiovascular functions. OX2 is one of the two receptors for the orexins. The receptor is expressed in the brain regions such as the hypothalamus, hippocampus and thalamus. Recent studies have indicated that orexins may also be useful in the treatment of schizophrenia and other psychiatric disorders.

Application: Functional assays

Figure 1

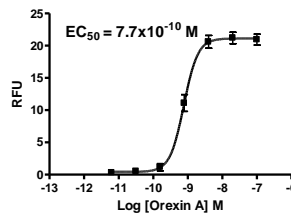


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Deutch and Bubser (2007) The orexins/hypocretins and schizophrenia. *Schizophr Bull* 33:1277-1283.

Smart *et al.* (1999) Characterization of recombinant human orexin receptor pharmacology in a Chinese hamster ovary cell-line using FLIPR. *Br J Pharmacol* 128:1-3.

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.