

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CXCR2 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** CG1002

**Lot Number:** CG1002-012313

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T Gaqi5

**Transfection:** Expression vector containing full-length human CXCR2 cDNA (GenBank Accession Number NM\_001557) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

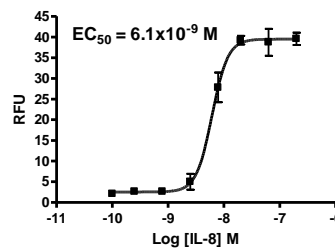
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin, 250  $\mu$ g/mL hygromycin

**Stability:** Testing in progress

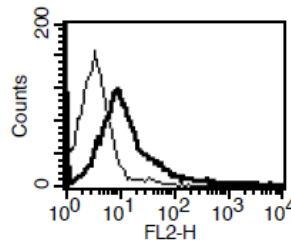
**Background:** CXCR2 (high affinity interleukin-8 receptor B, IL8RB) is a receptor to interleukin-8, which is a powerful neutrophil chemotactic factor. Binding of IL-8 to the receptor causes activation of neutrophils. This receptor also binds to GRO/MGSA and NAP-2 with high affinity. Some highly potent small molecule antagonists of CXCR2 showed success in blocking *in vivo* trafficking of neutrophils, suggesting that antagonism of IL-8 at the receptor level is a viable therapeutic strategy in conditions such as pulmonary disease.

**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:

Sprenger *et al.* (1994) Structure, genomic organization, and expression of the human interleukin-8 receptor B gene. *J Biol Chem* 269:11065-11072.

Pease and Sabroe (2002) The role of interleukin-8 and its receptors in inflammatory lung disease: implications for therapy. *Am J Respir Med* 1:19-25.

**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.