

## MULTISCREEN™ DIVISION ARRESTED CELL LINE HUMAN RECOMBINANT FPR2 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** DC1244a

**Lot Number:** DC1244a-121522

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cellbanker 2 (AmsBio 11891)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human FPR2 cDNA (GenBank Accession Number NM\_001462.3) with FLAG tag sequence at N-terminus

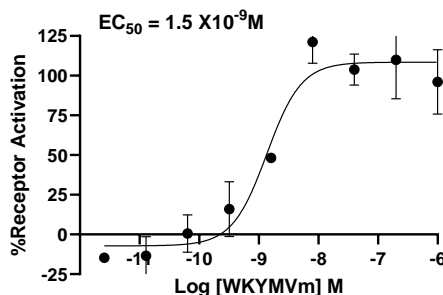
**Recommended Storage:** Liquid nitrogen upon receiving

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

**Background:** The lipoxin A4 receptor ALX is also known as formyl peptide receptor-like 1 (FPRL1) or formyl peptide receptor 2 (FPR2). It shares 69% amino acid identity with FPR but displays low affinity for bacterial peptide N-formyl-methionyl-leucyl-phenylalanine (fMLF). ALX is highly expressed on neutrophils and monocytes and mediates cell chemotaxis. By interacting with a variety of exogenous and host-derived agonists, ALX may have important implications in the pathogenesis of human diseases such as HIV, Alzheimer's disease (AD), amyloidosis and prion diseases.

**Application:** Functional assays

**Figure 1**



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

#### References:

Le *et al.* (2002) Formyl-peptide receptors revisited. *Trends Immunol* 23:541-548.

Migeotte *et al.* (2005) Identification and characterization of an endogenous chemotactic ligand specific for FPRL2. *J Exp Med* 201:83-93.

**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.