

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT C3A RECEPTOR**

PRODUCT INFORMATION

Catalog Number: DC1222

Lot Number: DC1222-090419

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

Freeze Medium: Cell Banker 2

Host cell: HEK293T

Transfection: Expression vector containing full-length human C3AR1 cDNA (GenBank Accession Number NM_004054.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Data sheet

Background: C3a receptor (C3a anaphylatoxin chemotactic receptor) is a receptor for the chemotactic and inflammatory peptide anaphylatoxin C3a. The receptor is widely expressed in many hematopoietic cell lines, throughout the CNS and other tissues, and the expression level in glial cells increases during CNS inflammation. It stimulates chemotaxis, granule enzyme release and superoxide anion production, mediates the pathogenesis of demyelinating disease by directly or indirectly chemoattracting encephalitogenic cells to the CNS. Administration of small molecule inhibitors of C3a receptor inhibits airway responses.

Application: Functional assays

Figure 1

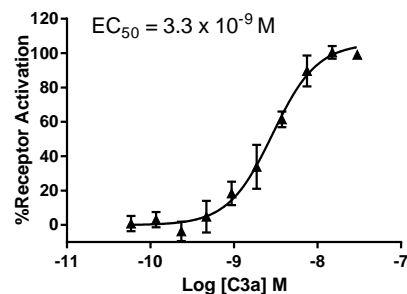


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Boos *et al.* (2004) Deletion of the complement anaphylatoxin C3a receptor attenuates, whereas ectopic expression of C3a in the brain exacerbates, experimental autoimmune encephalomyelitis. *J Immunol* 173:4708-14.

Ali and Panettieri, Jr. (2005) Anaphylatoxin C3a receptors in asthma. *Respir Res* 6:19.

FOR RESEARCH USE ONLY.

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.