

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT C5A RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1221

**Lot Number:** C1221-010106

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

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

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human C5aR1 cDNA (GenBank Accession Number NM\_001736) with FLAG tag sequence at N-terminus

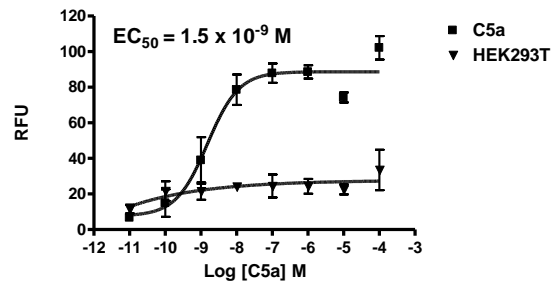
**Recommended Storage:** Liquid nitrogen upon receiving

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

**Stability:** In progress

**Background:** C5a receptor (C5a anaphylatoxin chemotactic receptor, or CD88 antigen, C5R1) is a receptor for chemotactic and inflammatory peptide anaphylatoxin C5a. C5a is a classic, potent inflammatory mediator, and the C5R1 receptor stimulates chemotaxis, granule enzyme release and superoxide anion production. C5a receptor antagonists, such as AcPhe[Orn-Pro-D-cyclohexylalanine-Trp-Arg], show therapeutic effects in various inflammatory disease models. C5a receptor may also be involved in neurodegeneration and Alzheimer's disease.

**Application:** Functional assay



**Figure legend:** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

#### References:

Sumichika, H. (2004). "C5a receptor antagonists for the treatment of inflammation." *Curr Opin Investig Drugs* 5(5): 505-10. Woodruff, T. M., T. V. Arumugam, et al. (2003). "A potent human C5a receptor antagonist protects against disease pathology in a rat model of inflammatory bowel disease." *J Immunol* 171(10): 5514-20.

Farkas, I., M. Takahashi, et al. (2003). "Complement C5a receptor-mediated signaling may be involved in neurodegeneration in Alzheimer's disease." *J Immunol* 170(11): 5764-71.

**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.