

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CRF1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** CG1040

**Lot Number:** CG1040-123013

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

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

**Host cell:** HEK293T Gαq5

**Transfection:** Expression vector containing full-length human CRHR1 cDNA (GenBank Accession Number NM\_004382.3) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

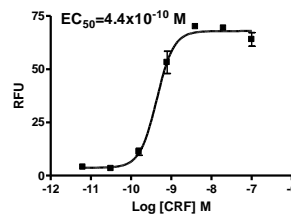
**Propagation Medium:** DMEM, 10% FBS, 1 μg/mL puromycin, 250 μg/mL hygromycin

**Stability:** Stable in culture for minimum of two months

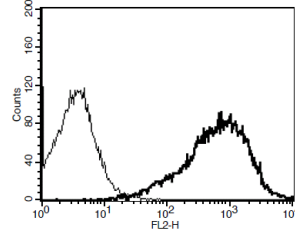
**Background:** Hypothalamic corticotropin releasing hormone (CRH) exerts its diverse physiological actions via cell surface G protein-coupled CRH receptors (CRHRs), of which two subtypes (CRHR1 and CRHR2) have been identified. CRH regulates pituitary ACTH secretion and mediates behavioral and autonomic responses to stress. Overproduction of CRH and stress system abnormalities are associated with psychiatric diseases such as depression, anxiety, eating disorders, and addiction. CRHR1 is expressed in pituitary corticotrophs and the brain. CRH binds to CRHR1 with high affinity, and activation of CRHR1 by CRH increases cAMP intracellular levels and activates protein kinase A. CRHR1 also couples to activation of MAPK as well as protein kinase C in an isoenzyme-specific manner. Selective CRHR1 antagonists offer new possibilities for the treatment of anxiety and depression.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

**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:

Dermitzaki *et al.* (2005) Corticotropin-releasing hormone activates protein kinase C in an isoenzyme-specific manner. *Biochem Biophys Res Commun* 327:828-836.

Muller and Wurst (2004) Getting closer to affective disorders: the role of CRH receptor systems. *Trends Mol Med* 10:409-415.

Refojo *et al.* (2005) Corticotropin-releasing hormone activates ERK1/2 MAPK in specific brain areas. *Proc Natl Acad Sci U S A* 102:6183-6188.

#### 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.