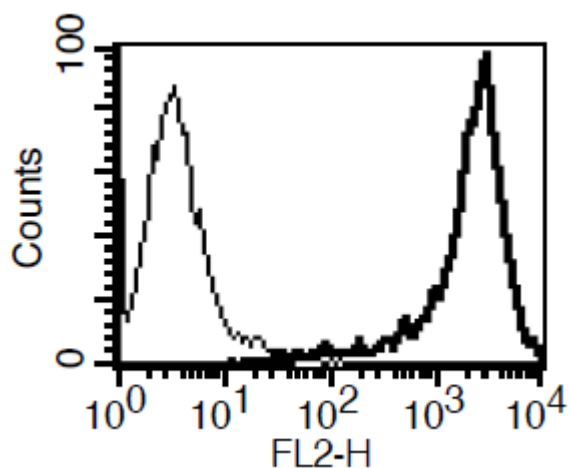


**MULTISCREEN™ STABLE CELL LINE MEMBRANE
RECOMBINANT HUMAN CASR RECEPTOR**

Materials Provided

Product Name: Human CASR Membrane	Cellular Background: HEK293T
Product Number: MC1233	Lot Number: MC1233-03202013
Protein Concentration 18.08 mg/mL	mL per vial 55 µL/vial
Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl ₂ , 25 mM Sucrose (pH 7.2)	
Storage Conditions: Store at -80°C. Avoid repeated freeze-thaw.	

Lot Specific Data



Histogram showing receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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. Mar 2010

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.