

MULTISCREEN™ HEK293T CELL LINE MEMBRANE

Data sheet

PRODUCT INFORMATION

Catalog Number: MC0001

Lot Number: MC0001-070215

Quantity: 17.3 mg/ml, 1 mg

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, 25 mM Sucrose (pH 7.2)

Host cell: HEK293T

Recommended Storage: Liquid nitrogen upon receiving

Figure 1:

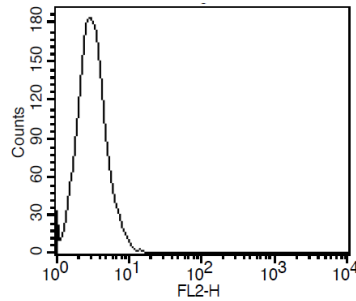


Figure 1: Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: HEK293T parental cells

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. October 2005

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