

**MULTISCREEN™ MEMBRANE PREPARATION
CHO-K1 PARENTAL CELLS**

Data sheet

PRODUCT INFORMATION

Catalog Number: MC0001-1

Lot Number: MC0001-1-052215

Quantity: 11.97 mg/ml, 1 mg

Packaging Buffer: 20 mM Gly-Gly,
1 mM MgCl₂, 250 mM Sucrose

Host cell: CHO-K1

Recommended Storage: Liquid
nitrogen upon receiving

Application: FACS

Figure 1

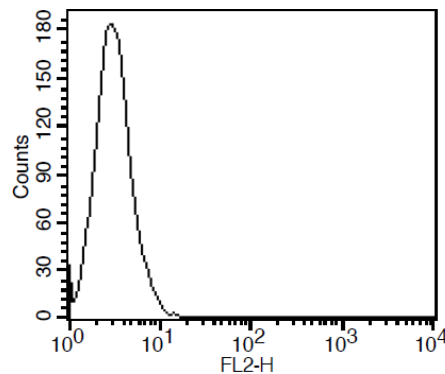


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells.

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.