

**MULTISCREEN™ MEMBRANE PREPARATION
HUMAN RH7777 PARENTAL**

Data sheet

PRODUCT INFORMATION

Catalog Number: MC0001-6

Lot Number: MC0001-6-042523

Quantity: 1 vial (9.7 mg/mL),
10mg or 1mg

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

Host cell: RH7777

Recommended Storage: Liquid
nitrogen upon receiving. Avoid
repeated freeze-thaw.

Figure 1.

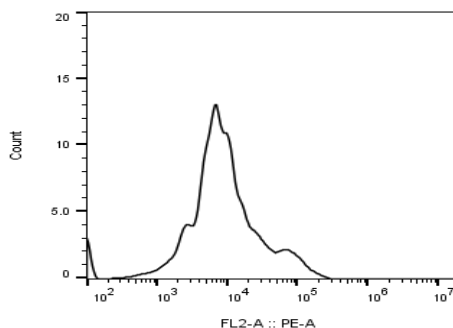


Figure 1: Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Black line: parental cells.

FOR RESEARCH USE ONLY.