

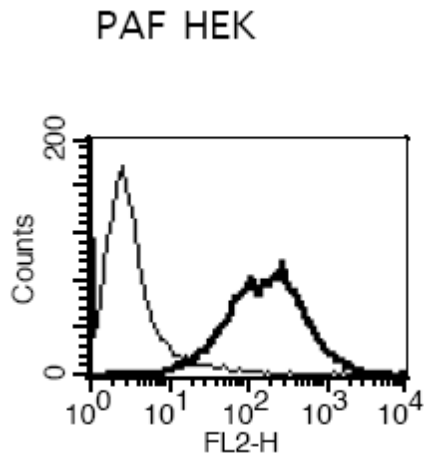
**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT PAF RECEPTOR MEMBRANE**

---

<b>Product Name:</b> PAF Membrane	<b>Cellular Background:</b> HEK293T
<b>Product Number:</b> MC1218	<b>Lot Number:</b> 03/25/2010
<b>Protein Concentration</b> 8.5 mg/mL	<b>Number of Vials/mL per vial</b> 12 vials, 1 mL/vial
<b>Packaging Buffer:</b> 20 mM Tris-HCl (pH7.5), 2 mM EDTA, protease inhibitor cocktail	
<b>Storage Conditions:</b> Store at -80°C. Avoid repeated freeze-thaw.	

**Lot Specific Data**

03/26/2010 FACS Analysis



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.