

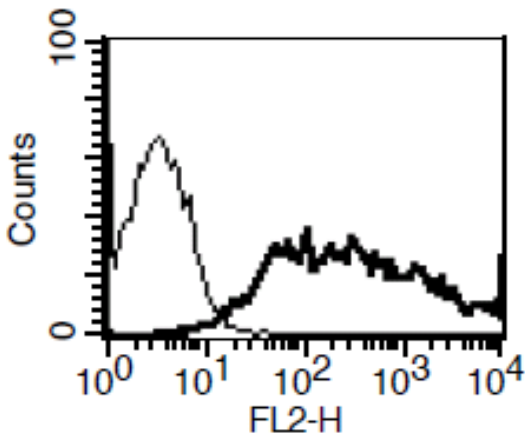
**MULTISCREEN™ STABLE CELL LINE MEMBRANE  
RECOMBINANT HUMAN ANGIOTENSIN II, TYPE 1 RECEPTOR**

---

**Materials Provided**

<b>Product Name:</b> AT1 Membrane	<b>Cellular Background:</b> HEK293T
<b>Product Number:</b> MH1417	<b>Lot Number:</b> MH1417-07152013
<b>Protein Concentration</b> 21.08 mg/mL	<b>mL per vial</b> 1 mg/vial
<b>Packaging Buffer:</b> 20 mM Gly-Gly, 1 mM MgCl <sub>2</sub> , 250 mM Sucrose	
<b>Storage Conditions:</b> 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.