

## MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT OT RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** MC1299-1

**Lot Number:** MC1299-1-111215

**Quantity:** 4.2mg/ml, 1mg

**Packaging Buffer:** 20 mM Gly-Gly, 1 mM MgCl<sub>2</sub>, 25 mM Sucrose (pH 7.2)

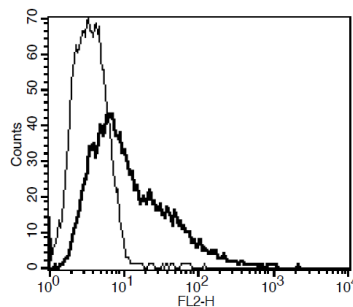
**Host cell:** CHO-K1

**Transfection:** Full-length Human OXTR cDNA (GenBank Accession Number NM\_000619.3) with FLAG-tag sequence at the N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving, avoid freeze-thaw

**Background:** Oxytocin (OT) is a member of the neurohypophyseal hormone family. OT receptors are found in uterine smooth muscle, myoepithelial cells in the mammary gland, and in the pituitary. OT stimulates contraction of uterine smooth muscle during labor and stimulates milk secretion in response to suckling. It has been postulated that OT may also facilitate social and bonding behaviors related to the reproduction and care of offspring. Further studies of the OT receptor are essential in gaining a better understanding of the mechanism and in vivo regulation of uterine function.

**Figure 1**



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

#### References:

Gimpl and Fahrenholz (2001) The oxytocin receptor system: structure, function, and regulation. *Physiol Rev* 81:629-683.

Ivell *et al.* (2001) The structure and regulation of the oxytocin receptor. *Exp Physiol* 86:289-296.

Shojo and Kaneko (2000) Characterization and expression of oxytocin and the oxytocin receptor. *Mol Genet Metab* 71:552-558.

**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.