

PRODUCT DATASHEET**ChemiScreen™ Wild-Type CHO Membrane Preparation**

CATALOG NUMBER:	HTS000MC	QUANTITY:	200 units
LOT NUMBER:	SC20190318	VOLUME/CONCENTRATION:	1 mL, 1 mg/mL
BACKGROUND:	CHO cells are commonly used for overexpression of GPCRs, and are used as a host for certain Chemicon ChemiScreen™ GPCR Membrane Preparations. Wild-Type CHO Membrane Preparations are ideal controls for validation of radioligand binding assays employing ChemiScreen™ GPCR Membrane Preparations in the CHO host cell line.		
APPLICATIONS:	Negative control in radioligand binding assay		
PRESENTATION:	Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA with no preservatives. Packaging method: Membrane proteins were adjusted to the indicated concentration in 1 ml packaging buffer, rapidly frozen, and stored at -80°C.		
STORAGE/HANDLING:	Store at -70°C. Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.		

FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

Eurofins Pharma Bioanalytics Services US Inc. is an independent member of Eurofins Discovery Services