

PRODUCT DATASHEET
ChemiScreen™ PAR2 Protease-Activated Membrane Preparation

CATALOG NUMBER: HTS037M **QUANTITY:** 200 units
LOT NUMBER: **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

BACKGROUND: The protease-activated receptor family of GPCRs has a unique mechanism of activation, in which protease cleaves a prodomain to reveal a peptide sequence that functions as a tethered ligand to activate the receptor (Macfarlane *et al.*, 2001). PAR2 is specifically activated by trypsin and mast cell tryptase, and can also be activated by free peptide analogs of the tethered ligand, such as SLIGRL and furoyl-LIGRLO (Coelho *et al.*, 2003; Kawabata *et al.*, 2004). PAR2 is expressed in endothelium, gastrointestinal epithelium, macrophages, eosinophils and nociceptive afferent neurons. Activation of PAR2 in these cells and tissues promotes vasodilation, inflammation, allergy, hyperalgesia and intestinal permeability. Therefore, PAR2 is regarded as an attractive therapeutic target for colitis, asthma, myocardial ischemia/reperfusion injury, and pain (Cocks *et al.*, 1999; Hansen *et al.*, 2005; Lindner *et al.*, 2000; McLean *et al.*, 2002; Vergnolle *et al.*, 2001). PAR2 membrane preparations are crude membrane preparations made from our proprietary stable recombinant cell lines to ensure high-level of GPCR surface expression; thus, they are ideal HTS tools for screening of antagonists of PAR2 interactions with its ligands. The membrane preparations exhibit an EC₅₀ of 1.9 μM for furoyl-LIGRLO in a GTP_γS binding assay.

APPLICATIONS: GTP_γS Binding

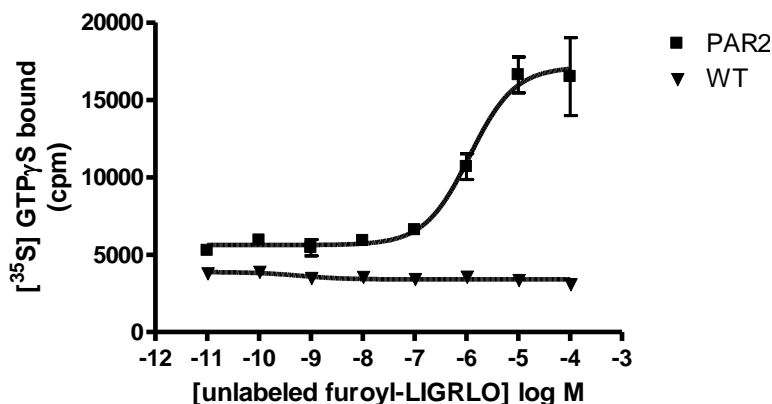


Figure 1. Binding of [³⁵S]GTP_γS to PAR2 membrane preparation. 5 μg/well PAR2 Membrane Preparation (catalog # HTS037M) and Wild-Type Chem-1 Membrane Preparation (catalog # HTS000MC1) was incubated with 0.1 nM [³⁵S]-GTP_γS and increasing amounts of unlabeled furoyl-LIGRLO. Bound radioactivity was determined by filtration and scintillation counting.

SPECIFICATIONS: 1 unit = 5 µg
EC50 for furoyl-LIGRLO: ~3.8 µM in a GTP γ S binding assay

Species: Human PAR2 (Accession number AY336105)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous PAR2 expression.

RECOMMENDED ASSAY CONDITIONS: Membranes are permeabilized by addition of saponin to an equal concentration by mass, then mixed with [³⁵S]-GTP γ S (final concentration of 0.1 nM) in 20 mM HEPES, pH 7.4/100 mM NaCl/10 mM MgCl₂/0.5 µM GDP in a nonbinding 96-well plate. Unlabeled furoyl-LIGRLO is added to the final concentration indicated in Figure 1 (final volume 100 µL), and incubated for 30 min at 30°C. The binding reaction is transferred to an FB filter plate (EMD Millipore MAHF B1H) prewetted with water, and washed 3 times (1 mL per well per wash) with cold 10 mM sodium phosphate, pH 7.4. The plate is dried and counted.

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than 1000 cpm specific furoyl-LIGRLO-stimulated [³⁵S]-GTP γ S binding.

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 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.

- REFERENCES:**
1. Cocks TM et al. (1999) A protective role for protease-activated receptors in the airways. *Nature* 398: 156-60.
 2. Coelho AM et al. (2003) Proteinase-activated receptor-2: physiological and pathophysiological roles. *Curr. Med. Chem. Cardiovasc. Hematol. Agents*. 1: 61-72.
 3. Hansen KK et al. (2005) A major role for proteolytic activity and proteinase-activated receptor-2 in the pathogenesis of infectious colitis. *Proc. Natl. Acad. Sci. USA*. 102: 8363-8.
 4. Kawabata et al. (2004) Potent and metabolically stable agonists for protease-activated receptor-2: evaluation of activity in multiple assay systems in vitro and in vivo. *J. Pharmacol. Exp. Ther.* 309: 1098-1107.
 5. Lindner JR et al. (2000) Delayed onset of inflammation in protease-activated receptor-2-deficient mice. *J. Immunol.* 165: 6504-10.
 6. Macfarlane SR et al. (2001) Proteinase-activated receptors. *Pharma. Rev.* 53: 245-82.
 7. McLean PG et al. (2002) Protease-activated receptor-2 activation causes EDHF-like coronary vasodilation: selective preservation in ischemia/reperfusion injury: involvement of lipoxigenase products, VR1 receptors, and C-fibers. *Circ. Res.* 90: 465-72.
 8. Vergnolle N et al. (2001) Proteinase-activated receptor-2 and hyperalgesia: A novel pain pathway. *Nat. Med.* 7: 821-6.

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