



MULTISPAN

Assay Experts in qPCRs and Beyond

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT EP4 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1204

Lot Number: DC1204-080625

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Cellbanker 2

Host cell: HEK293T

Transfection: Expression vector containing full-length human PTGER4 cDNA (GenBank Accession Number NM_000958) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Background: The human prostaglandin E2 (PGE2) receptor EP4 (or PTGER4) is abundantly expressed. Activation of EP4 receptor induces osteoblasts and thereby stimulates *de novo* bone formation. Results from targeted deletion of the EP4 receptor also suggest that the EP4 subtype of the PGE2 receptor is involved in lipopolysaccharide-induced bone resorption. Modulating EP4 action may thus prove to be clinically useful for the treatment of bacterially induced bone loss, such as in periodontitis and osteomyelitis. EP4 also plays a role in the anti-inflammatory action of PGE2 in human macrophage.

Application: Functional assays

Figure 1

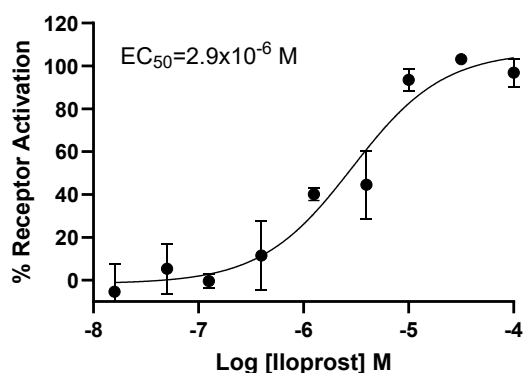


Figure 1. Dose-dependent accumulation of intracellular cAMP upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Morath *et al.* (1999) Immunolocalization of the four prostaglandin E2 receptor proteins EP1, EP2, EP3, and EP4 in human kidney. *J Am Soc Nephrol* 10:1851-1860.

Zhang *et al.* (2000) Characterization of murine vasopressor and vasodepressor prostaglandin E(2) receptors. *Hypertension* 35:1129-1134.

Li X *et al.* (2009) Prostaglandin E(2) and its cognate EP receptors control human adult articular cartilage homeostasis and are linked to the pathophysiology of osteoarthritis. *Arthritis Rheum* 60:513-523.

FOR RESEARCH USE ONLY.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

26219 Eden Landing Road, Hayward, California 94545-3718 Phone: (510) 887-0817 info@multispaninc.com