

# PRODUCT INFORMATION

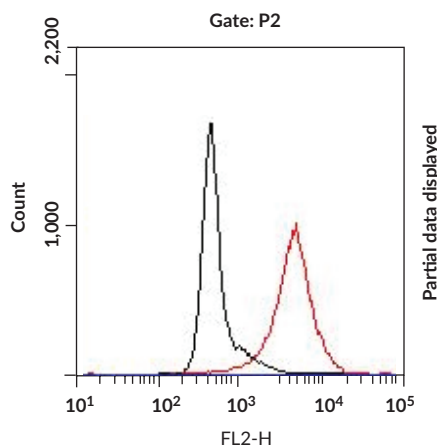


## EP<sub>2</sub> Receptor Polyclonal PE Antibody Item No. 10477

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of peptide affinity purified, PE labeled polyclonal antibody.
<b>Synonyms:</b>	PGE <sub>2</sub> Receptor 2, Prostaglandin E <sub>2</sub> Receptor 2
<b>Immunogen:</b>	Peptide from the C-terminal region of human EP <sub>2</sub>
<b>Cross Reactivity:</b>	(+) Human, mouse, and rat EP <sub>2</sub> receptors; (-) EP <sub>1</sub> , EP <sub>3</sub> , and EP <sub>4</sub> receptors
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Flow cytometry (FC) and immunocytochemistry (ICC); the recommended starting dilution for FC and ICC is 1:50. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Jurkat cell EP2 receptor detected with anti-EP2: R-phycoerythrin at 4 µg/ml

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/15/2024

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

This polyclonal antibody is labeled with R-Phycoerythrin (PE) and can be used for immunofluorescent labeling of cellular EP<sub>2</sub> receptor following permeabilization. R-PE absorption maxima occurs at 565>540>498 nm and the emission maximum occurs at 578 nm, however standard excitation by 488 nm lasers is sufficient to generate signal detected in the PE channel of flow cytometers. The biological effects of prostaglandin E<sub>2</sub> (PGE<sub>2</sub>) are mediated through interaction with four distinct membrane-bound G-protein coupled EP receptors: EP<sub>1</sub>, EP<sub>2</sub>, EP<sub>3</sub>, and EP<sub>4</sub>.<sup>1,2</sup> Binding of PGE<sub>2</sub> to the EP<sub>2</sub> receptor results in an increase in adenylate cyclase activity with a subsequent increase in cAMP.<sup>3,4</sup> Pharmacologically, EP<sub>2</sub> receptors mediate relaxation of smooth muscle and are distinguished from EP<sub>4</sub> receptors by their sensitivity to the EP<sub>2</sub> receptor selective agonist butaprost.<sup>1-3</sup> The human EP<sub>2</sub> receptor is comprised of 358 amino acids with a molecular mass of approximately 40,000.<sup>3</sup> mRNA for the EP<sub>2</sub> receptor is expressed in a variety of tissues including lung, placenta, spleen, intestine, kidney, and sensory neuron.<sup>3-5</sup>

## References

---

1. Coleman, R.A., Smith, W.L., and Narumiya, S. International Union of Pharmacology classification of prostanoid receptors: Properties, distribution, and structure of the receptors and their subtypes. *Pharmacol. Rev.* **46(2)**, 205-229 (1994).
2. Coleman, R.A., Eglén, R.M., Jones, R.L., *et al.* Prostanoid and leukotriene receptors: A progress report from the IUPHAR working parties on classification and nomenclature. *IUPHAR Compendium* 1-12 (1997).
3. Narko, K., Saukkonen, K., Ketola, I., *et al.* Regulated expression of prostaglandin E<sub>2</sub> receptors EP<sub>2</sub> and EP<sub>4</sub> in human ovarian granulosa-luteal cells. *J. Clin. Endocrinol. Metab.* **86(4)**, 1765-1768 (2001).
4. Nemoto, K., Pilbeam, C.C., Bilak, S.R., *et al.* Molecular cloning and expression of a rat prostaglandin E<sub>2</sub> receptor of the EP<sub>2</sub> subtype. *Prostaglandins* **54(4)**, 713-725 (1997).
5. Southall, M.D. and Vasko, M.R. Prostaglandin receptor subtypes, EP<sub>3C</sub> and EP<sub>4</sub>, mediate the prostaglandin E<sub>2</sub>-induced cAMP production and sensitization of sensory neurons. *J. Biol. Chem.* **276(19)**, 16083-16091 (2001).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM