

# PRODUCT INFORMATION

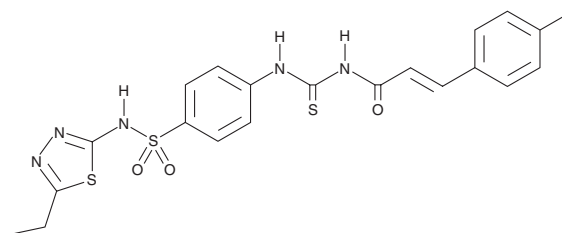


## TG6-129

Item No. 19085

**CAS Registry No.:** 1164464-14-5  
**Formal Name:** (2E)-N-[[[4-[[[(5-ethyl-1,3,4-thiadiazol-2-yl)amino]sulfonyl]phenyl]amino]thioxomethyl]-3-(4-fluorophenyl)-2-propenamide

**Synonym:** SID 17503974  
**MF:** C<sub>20</sub>H<sub>18</sub>FN<sub>5</sub>O<sub>3</sub>S<sub>3</sub>  
**FW:** 491.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 310 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

TG6-129 is supplied as a crystalline solid. A stock solution may be made by dissolving the TG6-129 in the solvent of choice, which should be purged with an inert gas. TG6-129 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TG6-129 in these solvents is approximately 10 and 8 mg/ml, respectively.

TG6-129 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010) evokes distinct responses through four different 'E prostanoid' (EP) receptors. EP<sub>2</sub> is a G protein-coupled receptor that has diverse roles, including those in cancer, inflammation, and neuroprotection.<sup>1-3</sup> TG6-129 is an antagonist of the EP<sub>2</sub> receptor, suppressing PGE<sub>2</sub>-induced elevation of cAMP in cells expressing EP<sub>2</sub> with an IC<sub>50</sub> value of 1.6 μM.<sup>4</sup> It is without effect on EP<sub>4</sub>, DP<sub>1</sub>, IP, and β<sub>2</sub>-adrenergic receptors. TG6-129 reduces the expression of COX-2, IL-1β, IL-12, IL-23, IL-6, and TNF-α induced by the EP<sub>2</sub>-selective agonist butaprost (Item No. 13740) in P388D1 macrophages.<sup>4</sup> It has low cell cytotoxicity (CC<sub>50</sub> = 326 μM), prolonged plasma half-life, and does not cross the blood-brain barrier.<sup>4</sup>

### References

1. Majima, M., Amano, H., and Hayashi, I. Prostanoid receptor signaling relevant to tumor growth and angiogenesis. *Trends Pharmacol. Sci.* **34**(10), 524-529 (2003).
2. Jiang, J. and Dingleline, R. Prostaglandin receptor EP<sub>2</sub> in the crosshairs of anti-inflammation, anti-cancer, and neuroprotection. *Trends Pharmacol. Sci.* **34**(7), 413-423 (2013).
3. Kawahara, K., Hohjoh, H., Inazumi, T., et al. Prostaglandin E<sub>2</sub>-induced inflammation: Relevance of prostaglandin E receptors. *Biochim. Biophys. Acta* **1851**(4), 414-421 (2015).
4. Ganesh, T., Jiang, J., Shashidharamurthy, R., et al. Discovery and characterization of carbamothioylacrylamides as EP<sub>2</sub> selective antagonists. *ACS Med. Chem. Lett.* **4**(7), 616-621 (2013).

#### 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, 11/28/2022

#### 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