

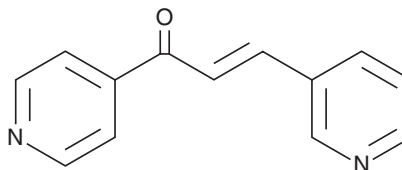
# PRODUCT INFORMATION



**3PO**

Item No. 19276

**CAS Registry No.:** 18550-98-6  
**Formal Name:** 3-(3-pyridinyl)-1-(4-pyridinyl)-2E-propen-1-one  
**MF:** C<sub>13</sub>H<sub>10</sub>N<sub>2</sub>O  
**FW:** 210.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 226, 302 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

3PO is supplied as a crystalline solid. A stock solution may be made by dissolving the 3PO in the solvent of choice, which should be purged with an inert gas. 3PO is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 3PO in these solvents is approximately 2, 25, and 5 mg/ml, respectively.

3PO is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3PO should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 3PO has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

3PO is a competitive inhibitor of 6-phosphofructo-2-kinase/fructose-2,6-bisphosphate 3 (PFKFB3; IC<sub>50</sub> = 23 μM) that causes a rapid reduction in fructose-2,6-bisphosphate (F2,6BP), glucose uptake, and lactate secretion.<sup>1,2</sup> This is followed by a reduction in the steady-state concentration of ATP and NADH, resulting in cell cycle arrest in Jurkat T cell leukemia cells.<sup>1,2</sup> 3PO is selectively cytostatic to *ras*-transformed bronchial epithelial cells relative to normal bronchial epithelial cells.<sup>1</sup> 3PO markedly reduces intracellular F2,6BP, glucose uptake, and growth of established tumors in mice.<sup>1</sup> It also blocks angiogenesis, reducing vessel sprouting in endothelial cell spheroids, zebrafish embryos, and the postnatal mouse retina by inhibiting endothelial cell proliferation and migration.<sup>3</sup>

## References

1. Clem, B., Telang, S., Clem, A., *et al.* Small-molecule inhibition of 6-phosphofructo-2-kinase activity suppresses glycolytic flux and tumor growth. *Mol. Cancer Ther.* **7**(1), 110-120 (2008).
2. Clem, B.F., O'Neal, J., Tapolsky, G., *et al.* Targeting 6-phosphofructo-2-kinase (PFKFB3) as a therapeutic strategy against cancer. *Mol. Cancer Ther.* **12**(8), 1461-1470 (2013).
3. Schoors, S., De Bock, K., Cantelmo, A.R., *et al.* Partial and transient reduction of glycolysis by PFKFB3 blockade reduces pathological angiogenesis. *Cell Metab.* **19**(1), 37-48 (2014).

### 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, 12/01/2022

## 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](http://WWW.CAYMANCHEM.COM)