

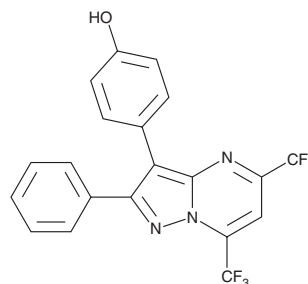
# PRODUCT INFORMATION



## PHTPP

Item No. 16025

**CAS Registry No.:** 805239-56-9  
**Formal Name:** 4-[2-phenyl-5,7-bis(trifluoromethyl)pyrazolo[1,5-a]pyrimidin-3-yl]-phenol  
**MF:** C<sub>20</sub>H<sub>11</sub>F<sub>6</sub>N<sub>3</sub>O  
**FW:** 423.3  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

PHTPP is supplied as a crystalline solid. A stock solution may be made by dissolving the PHTPP in the solvent of choice, which should be purged with an inert gas. PHTPP is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PHTPP in ethanol is approximately 30 mg/ml and approximately 50 mg/ml in DMSO and DMF.

PHTPP is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PHTPP should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PHTPP has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

PHTPP is a pyrazolo[1,5- $\alpha$ ]pyrimidine-based ligand that acts as a full antagonist of estrogen ER $\beta$  receptors with 36-fold selectivity over ER $\alpha$ .<sup>1</sup> It exhibits no significant agonism on ER $\alpha$  or ER $\beta$ .<sup>1</sup> This compound has been used to selectively target ER $\beta$  in the study of the opposing effects of hormone therapy on tumors expressing either ER subtype. At 100 pM, PHTPP has been found to enhance SKOV3 and OV2008 ovarian cancer cell growth in *in vitro* assays.<sup>2</sup>

### References

1. Compton, D.R., Sheng, S., Carlson, K.E., *et al.* Pyrazolo[1,5- $\alpha$ ]pyrimidines: Estrogen receptor ligands possessing estrogen receptor  $\beta$  antagonist activity. *J. Med. Chem.* **47**(24), 5872-5893 (2004).
2. Chan, K.K.L., Leung, T.H.Y., Chan, D.W., *et al.* Targeting estrogen receptor subtypes (ER $\alpha$  and ER $\beta$ ) with selective ER modulators in ovarian cancer. *J. Endocrinol.* **221**(2), 325-336 (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, 02/14/2022

#### 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