

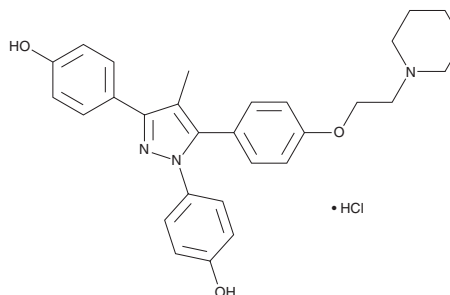
# PRODUCT INFORMATION



## Methylpiperidino pyrazole (hydrochloride)

Item No. 45218

**CAS Registry No.:** 2863676-89-3  
**Formal Name:** 4-[1-(4-hydroxyphenyl)-4-methyl-5-[4-[2-(1-piperidinyl)ethoxy]phenyl]-1H-pyrazol-3-yl]-phenol, monohydrochloride  
**Synonym:** MPP  
**MF:** C<sub>29</sub>H<sub>31</sub>N<sub>3</sub>O<sub>3</sub> • HCl  
**FW:** 506.0  
**Purity:** ≥95%  
**Supplied as:** A 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

Methylpiperidino pyrazole (MPP) (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the MPP (hydrochloride) in the solvent of choice, which should be purged with an inert gas. MPP (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

MPP is a selective estrogen receptor modulator (SERM).<sup>1,2</sup> It is an estrogen receptor  $\alpha$  (ER $\alpha$ ) antagonist ( $K_i = 2.7$  nM) and is selective for ER $\alpha$  over ER $\beta$  ( $K_i = 1,800$  nM).<sup>1</sup> MPP (1 nM) induces apoptosis in RL-95 endometrial cancer cells.<sup>2</sup> *In vivo*, MPP (25, 50, and 100  $\mu$ g/animal) increases uterine weight in ovariectomized mice, indicating differential *in vitro* versus *in vivo* endometrial responses.

### References

1. Sun, J., Huang, Y.R., Harrington, W.R., *et al.* Antagonists selective for estrogen receptor  $\alpha$ . *Endocrinology* **143**(3), 941-947 (2002).
2. Davis, A.M., Ellersieck, M.R., Grimm, K.M., *et al.* The effects of the selective estrogen receptor modulators, methyl-piperidino-pyrazole (MPP), and raloxifene in normal and cancerous endometrial cell lines and in the murine uterus. *Mol. Reprod. Dev.* **73**(8), 1034-1044 (2006).

#### 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, 03/03/2026

#### 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