

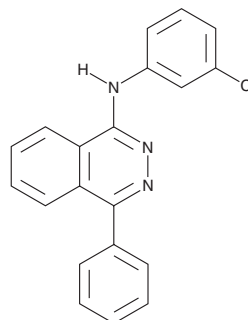
# PRODUCT INFORMATION



**MY-5445**

Item No. 20872

**CAS Registry No.:** 78351-75-4  
**Formal Name:** N-(3-chlorophenyl)-4-phenyl-1-phthalazinamine  
**MF:** C<sub>20</sub>H<sub>14</sub>ClN<sub>3</sub>  
**FW:** 331.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 215, 336 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

MY-5445 is supplied as a crystalline solid. A stock solution may be made by dissolving the MY-5445 in the solvent of choice, which should be purged with an inert gas. MY-5445 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of MY-5445 in these solvents is approximately 2 mg/ml.

MY-5445 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MY-5445 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MY-5445 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

MY-5445 is a selective inhibitor of the cyclic GMP (cGMP) phosphodiesterase, PDE5, with a K<sub>i</sub> value of 1.3 μM.<sup>1,2</sup> It inhibits calcium- and cAMP-selective PDEs with K<sub>i</sub> values of >1,000 and 915 μM, respectively.<sup>1</sup> MY-5445 has been shown to relax intact strips of rat aorta with an EC<sub>50</sub> value of 4 μM.<sup>1</sup>

## References

1. Souness, J.E., Brazdil, R., Diocee, B.K., *et al.* Role of selective cyclic GMP phosphodiesterase inhibition in the myorelaxant actions of M&B 22,948, MY-5445, vinpocetine and 1-methyl-3-isobutyl-8-(methylamino)xanthine. *Br. J. Pharmacol.* **98(3)**, 725-734 (1989).
2. Yamamoto, T., Yamamaoto, S., Osborne, Jr., J.C., *et al.* Complex effects of inhibitors on cyclic GMP-stimulated cyclic nucleotide phosphodiesterase. *J. Biol. Chem.* **258(23)**, 14173-14177 (1983).

### 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, 12/06/2022

## 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