

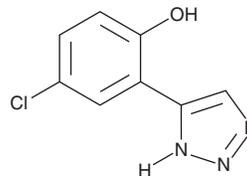
# PRODUCT INFORMATION



## MMG-0358

Item No. 29828

**CAS Registry No.:** 1378976-02-3  
**Formal Name:** 4-chloro-2-(1H-1,2,3-triazol-5-yl)-phenol  
**MF:** C<sub>8</sub>H<sub>6</sub>ClN<sub>3</sub>O  
**FW:** 195.6  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 214, 247, 304 nm  
**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

MMG-0358 is supplied as a solid. A stock solution may be made by dissolving the MMG-0358 in the solvent of choice, which should be purged with an inert gas. MMG-0358 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MMG-0358 in ethanol and DMSO is approximately 15 mg/ml and approximately 20 mg/ml in DMF.

MMG-0358 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MMG-0358 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MMG-0358 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

MMG-0358 is an inhibitor of indoleamine 2,3-dioxygenase 1 (IDO1; IC<sub>50</sub>s = 2 and 80 nM for the mouse and human enzyme, respectively).<sup>1,2</sup> It is selective for IDO1 over tryptophan 2,3-dioxygenase (TDO; IC<sub>50</sub>s = >100 μM for mouse and human TDO).<sup>1</sup>

### References

1. Röhrig, U.F., Majjigapu, S.R., Grosdidier, A., *et al.* Rational design of 4-aryl-1,2,3-triazoles for indoleamine 2,3-dioxygenase 1 inhibition. *J. Med. Chem.* **55**(11), 5270-5290 (2012).
2. Röhrig, U.F., Reynaud, A., Majjigapu, S.R., *et al.* Inhibition mechanisms of indoleamine 2,3-dioxygenase 1 (IDO1). *J. Med. Chem.* **62**(19), 8784-8795 (2019).

#### 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, 12/08/2022

#### 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