

# PRODUCT INFORMATION

## 1400W (hydrochloride)

Item No. 81520



CAS Registry No.: 214358-33-5

Formal Name: N-[[3-(aminomethyl)phenyl]methyl]-ethanimidamide, dihydrochloride

MF:  $C_{10}H_{15}N_3 \cdot 2HCl$

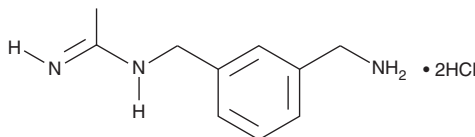
FW: 250.2

Purity:  $\geq 98\%$

Supplied as: A crystalline solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

1400W (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1400W (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 1400W (hydrochloride) is soluble in methanol and DMSO.

### Description

1400W is a potent, selective inhibitor of iNOS.<sup>1</sup> Using purified human iNOS, nNOS, and eNOS, 1400W exhibits  $K_i$  values of 7 nM, 2  $\mu$ M, and 50  $\mu$ M, respectively.<sup>1</sup> In rat aortic rings it is at least 1,000-fold more potent against iNOS compared to eNOS.<sup>1</sup> In mice bearing a murine mammary carcinoma, continuous infusion of 1400W at a rate of 10 mg/kg for 6 days decreased tumor weight approximately 50% compared to controls.<sup>2</sup>

### References

1. Garvey, E.P., Oplinger, J.A., Furfine, E.S., *et al.* 1400W is a slow, tight binding, and highly selective inhibitor of inducible nitric-oxide synthase *in vitro* and *in vivo*. *J. Biol. Chem.* **272**(8), 4959-4963 (1997).
2. Thomsen, L.L., Scott, J.M.J., Topley, P., *et al.* Selective inhibition of inducible nitric oxide synthase inhibits tumor growth *in vivo*: Studies with 1400W, a novel inhibitor. *Cancer Res.* **57**(15), 3300-3304 (1997).

#### 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, 06/29/2023

#### 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