

# PRODUCT INFORMATION



## CWHM12

Item No. 19480

**CAS Registry No.:** 1564286-55-0  
**Formal Name:** (3S)-N-[3-hydroxy-5-[(1,4,5,6-tetrahydro-5-hydroxy-2-pyrimidinyl)amino]benzoyl]glycyl-3-[3-bromo-5-(1,1-dimethylethyl)phenyl]-β-alanine

**MF:** C<sub>26</sub>H<sub>32</sub>BrN<sub>5</sub>O<sub>6</sub>

**FW:** 590.5

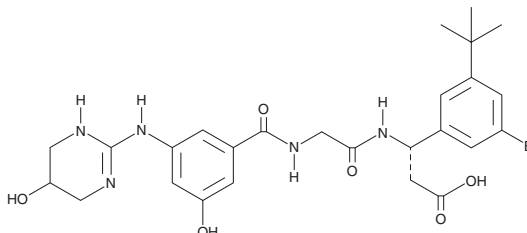
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 299 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

CWHM12 is supplied as a crystalline solid. A stock solution may be made by dissolving the CWHM12 in the solvent of choice. CWHM12 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of CWHM12 in these solvents is approximately 1 mg/ml.

CWHM12 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CWHM12 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CWHM12 has a solubility of approximately 0.33 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

CWHM12 is an analog of RGD peptide (Item No. 14501), a tripeptide that inhibits integrin-ligand interactions in studies related to cell adhesion, migration, growth, and differentiation.<sup>1</sup> CWHM12 selectively inhibits αv integrins (IC<sub>50</sub>s = 1.8, 0.8, 61, 1.5, and 0.2 nM for αvβ1, αvβ3, αvβ5, αvβ6, and αvβ8, respectively) over α11β3, α2β1, and α10β1.<sup>2</sup> CWHM 12 attenuates liver, lung, and pancreatic fibrosis in mice treated with CCl<sub>4</sub> or cerulein.<sup>2,3</sup>

### References

1. Ruoslahti, E. and Öbrink, B. Common principles in cell adhesion. *Exp. Cell Res.* **227**, 1-11 (1996).
2. Henderson, N.C., Arnold, T.D., Katamura, Y., et al. Selective αv integrin depletion identifies a core, targetable molecular pathway that regulates fibrosis across solid organs. *Nat. Med.* **19(12)**, 1-12 (2013).
3. Ulmasov, B., Newschwander-Tetri, B.A., Lai, J., et al. Inhibitors of Arg-Gly-Asp-binding integrins reduce development of pancreatic fibrosis in mice. *Cell. Mol. Gastroenterol. Hepatol.* **2(4)**, 499-518 (2016).

#### 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, 10/31/2022

#### 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