

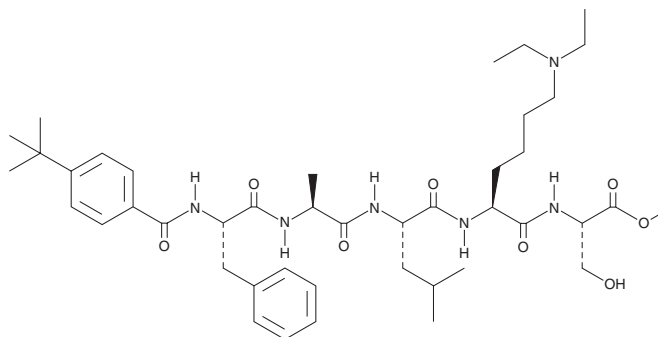
# PRODUCT INFORMATION



## UNC3866

Item No. 19237

**CAS Registry No.:** 1872382-47-2  
**Formal Name:** N-[4-(1,1-dimethylethyl)benzoyl]-L-phenylalanyl-L-alanyl-L-leucyl-N<sup>6</sup>,N<sup>6</sup>-diethyl-L-lysyl-L-serine, methyl ester  
**MF:** C<sub>43</sub>H<sub>66</sub>N<sub>6</sub>O<sub>8</sub>  
**FW:** 795.0  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 240 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

UNC3866 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC3866 in the solvent of choice. UNC3866 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of UNC3866 in ethanol is approximately 16 mg/ml and approximately 33 mg/ml in DMSO and DMF.

UNC3866 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNC3866 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. UNC3866 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

UNC3866 is an inhibitor of chromobox homolog 7 (CBX7; K<sub>d</sub> = 97 nM).<sup>1</sup> It also binds to CBX4 (K<sub>d</sub> = 94 nM) but is greater than 6-fold selective over CBX2, 6, and 8 and the chromodomain Y family (CDY) chromodomains CDY1, L1b, and L2 (K<sub>d</sub>s = 0.61-6.3 μM). UNC3866 pretreatment (30 μM) of PC3 lysates completely inhibits chemiprecipitation of CBX7. It induces a senescent-like morphology and reduces growth of PC3 cells with an EC<sub>50</sub> value of 7.6 μM.

### Reference

1. Stuckey, J.I., Dickson, B.M., Cheng, N., *et al.* A cellular chemical probe targeting the chromodomains of Polycomb repressive complex 1. *Nat. Chem. Biol.* **12**(3), 180-187 (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, 12/16/2022

#### 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