

# PRODUCT INFORMATION



**PD 173212**

Item No. 29447

**CAS Registry No.:** 217171-01-2  
**Formal Name:** N-[[4-(1,1-dimethylethyl)phenyl]methyl]-N-methyl-L-leucyl-N-(1,1-dimethylethyl)-O-(phenylmethyl)-L-tyrosinamide

**MF:** C<sub>38</sub>H<sub>53</sub>N<sub>3</sub>O<sub>3</sub>

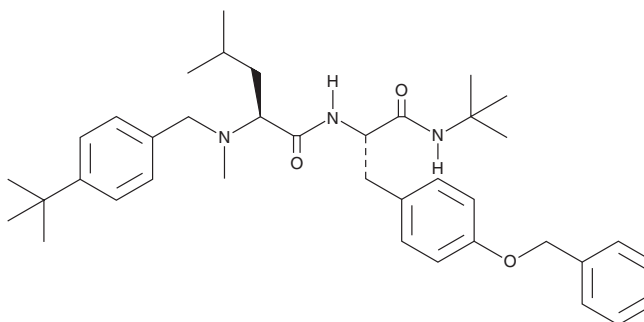
**FW:** 599.9

**Purity:** ≥95%

**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

PD 173212 is supplied as a solid. A stock solution may be made by dissolving the PD 173212 in the solvent of choice, which should be purged with an inert gas. PD 173212 is soluble in organic solvents such as ethanol and DMSO. The solubility of PD 173212 in these solvents is approximately 100 mM.

## Description

PD 173212 is an N-type calcium channel blocker (IC<sub>50</sub> = 36 nM).<sup>1</sup> It prevents audiogenic seizures in mice when administered at a dose of 30 mg/kg.

## Reference

1. Hu, L.-Y., Ryder, T.R., Rafferty, M.F., *et al.* Structure-activity relationship of N-methyl-N-aryl-peptidylamines as novel N-type calcium channel blockers. *Bioorg. Med. Chem. Lett.* **9(15)**, 2151-2156 (1999).

### 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, 01/03/2024

## 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](http://WWW.CAYMANCHEM.COM)