

# PRODUCT INFORMATION



## PKCε Inhibitor Scramble Peptide

Item No. 31761

**CAS Registry No.:** 813416-34-1  
**Formal Name:** L-leucyl-L-seryl-L-α-glutamyl-L-threonyl-L-lysyl-L-prolyl-L-alanyl-L-valine  
**Synonyms:** H-Leu-Ser-Glu-Thr-Lys-Pro-Ala-Val-OH, H-LSETKPAV-OH  
**MF:** C<sub>37</sub>H<sub>65</sub>N<sub>9</sub>O<sub>13</sub> H—Leu—Ser—Glu—Thr—Lys—Pro—Ala—Val—OH  
**FW:** 844.0  
**Purity:** ≥95%  
**Supplied as:** A lyophilized solid  
**Storage:** -20°C  
**Stability:** ≥3 years

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

### Laboratory Procedures

PKCε inhibitor scramble peptide is supplied as a lyophilized solid. A stock solution may be made by dissolving the PKCε inhibitor scramble peptide in water. The solubility of PKCε inhibitor scramble peptide in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PKCε inhibitor scramble peptide is a scrambled peptide with identical amino acid composition to PKCε inhibitor peptide (Item No. 17476). It is intended for use as a negative control for PKCε inhibitor peptide activity.

#### 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, 12/13/2023

#### 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