

PRODUCT INFORMATION



Acetyl Pentapeptide-1

Item No. 27261

CAS Registry No.:	97530-32-0
Formal Name:	N ² -acetyl-L-arginyl-L-lysyl-L- α -aspartyl-L-valyl-L-tyrosine
MF:	C ₃₂ H ₅₁ N ₉ O ₁₀
FW:	721.8
Purity:	≥95%
UV/Vis.:	λ_{max} : 276 nm
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

Acetyl pentapeptide-1 is supplied as a solid. Aqueous solutions of acetyl pentapeptide-1 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of acetyl pentapeptide-1 in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Acetyl pentapeptide-1 is a pentapeptide that decreases IL-8 secretion in human keratinocytes when used in combination with acetyl hexapeptide-36 and acetyl hexapeptide-38.¹ Formulations containing acetyl pentapeptide-1 have been used to decrease skin irritation caused by cleansing and as moisturizing agents in cosmetic applications.

Reference

1. Tittl, J. and Wolf, C. Compositions and methods for mitigating skin irritation and enhancing skin barrier function. *Gojo Industries, Inc. US 2016/0000858 A1* (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, 11/30/2022

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