

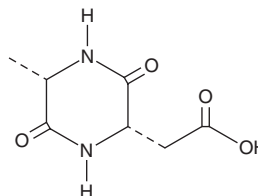
PRODUCT INFORMATION



Aspartyl-alanyl-diketopiperazine

Item No. 37715

CAS Registry No.: 110954-19-3
Formal Name: 5S-methyl-3,6-dioxo-2S-piperazineacetic acid
Synonym: DA-DKP
MF: C₇H₁₀N₂O₄
FW: 186.2
Purity: ≥90%
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

Aspartyl-alanyl-diketopiperazine is supplied as a solid. A stock solution may be made by dissolving the aspartyl-alanyl-diketopiperazine in the solvent of choice, which should be purged with an inert gas. Aspartyl-alanyl-diketopiperazine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of aspartyl-alanyl-diketopiperazine in these solvents is approximately 30 mg/ml. Aspartyl-alanyl-diketopiperazine is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of aspartyl-alanyl-diketopiperazine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of aspartyl-alanyl-diketopiperazine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Aspartyl-alanyl-diketopiperazine is an immunosuppressive diketopiperazine derivative of a human serum albumin (HSA) peptide.¹ It reduces antigen-induced TNF- α and IFN- γ production in isolated human T cells when used at a concentration of 500 μ M.

Reference

1. Shimonkevitz, R., Thomas, G., Slone, D.S., *et al.* A diketopiperazine fragment of human serum albumin modulates T-lymphocyte cytokine production through Rap1. *J. Trauma* **64**(1), 35-41 (2008).

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, 12/21/2023

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