

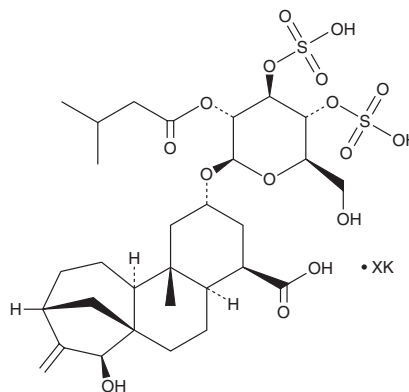
PRODUCT INFORMATION



Atractyloside (potassium salt)

Item No. 44522

CAS Registry No.: 695179-73-8
Formal Name: (2 β)-15 α -hydroxy-2-[[2-O-(3-methyl-1-oxobutyl)-3,4-di-O-sulfo- β -D-glucopyranosyl]oxy]-19-norkaur-16-en-18-oic acid, potassium salt
MF: C₃₀H₄₆O₁₆S₂ • XK
FW: 726.8
Purity: $\geq$ 95%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Laboratory Procedures

Atractyloside (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the atractyloside (potassium salt) in the solvent of choice, which should be purged with an inert gas. Atractyloside (potassium salt) is slightly soluble (0.1-1 mg/ml) in methanol and DMSO.

Description

Atractyloside is a natural heteroglucoside produced in some plants, including *A. gummifera*.¹ Atractyloside prevents mitochondrial ATP synthesis by inhibiting ADP/ATP translocases, which are responsible for the exchange of adenine di- and triphosphates across the inner mitochondrial membrane.^{1,2}

References

1. Dahout-Gonzalez, C., Hury, H., Trézéguet, V., *et al.* Molecular, functional, and pathological aspects of the mitochondrial ADP/ATP carrier. *Physiology (Bethesda)* **21**, 242-249 (2006).
2. Klingenberg, M. The ADP and ATP transport in mitochondria and its carrier. *Biochim. Biophys. Acta* **1778(10)**, 1978-2021 (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, 11/03/2025

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