

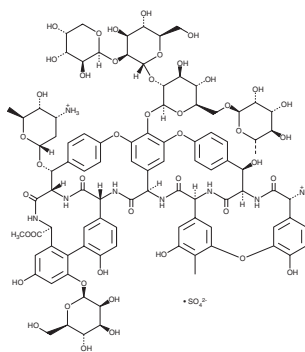
PRODUCT INFORMATION



Ristocetin A (sulfate)

Item No. 20601

CAS Registry No.: 90831-71-3
Formal Name: ristomycin A, monosulfate
Synonym: Ristomycin III
MF: $C_{95}H_{112}N_8O_{44} \cdot SO_4$
FW: 2,166.0
Purity: $\geq 95\%$
Supplied as: A powder
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Ristocetin A (sulfate) is supplied as a powder. A stock solution may be made by dissolving the ristocetin A (sulfate) in the solvent of choice, which should be purged with an inert gas. Ristocetin A (sulfate) is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Ristocetin A is an antibiotic that was used to treat staphylococcal infections.¹ However, side effects of ristocetin A included thrombocytopenia and platelet agglutination. Ristocetin A is now used in two assays related to these actions, the ristocetin cofactor assay and the ristocetin-induced platelet aggregation assay. Both assays are used in the diagnosis of von Willebrand disease and other bleeding disorders.^{1,2}

References

1. Hawks, G.H. Antibiotic therapy of staphylococcal infection. *Can. Med. Assoc. J.* **93**(16), 848-853 (1965).
2. Castaman, G., Hillarp, A., and Goodeve, A. Laboratory aspects of von Willebrand disease: Test repertoire and options for activity assays and genetic analysis. *Haemophilia* **20**(suppl 4), 65-70 (2014).

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, 10/04/2022

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