

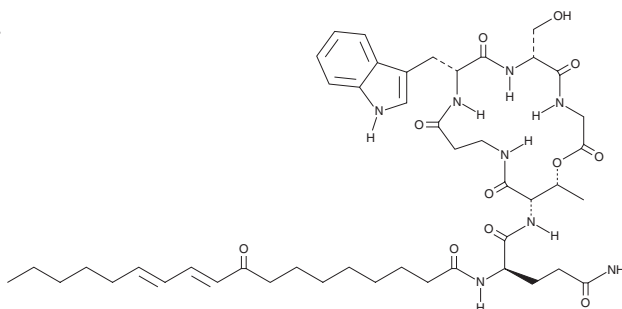
# PRODUCT INFORMATION



## Aselacin C

Item No. 33163

**CAS Registry No.:** 156223-08-4  
**Formal Name:** N<sup>2</sup>-(1,9-dioxo-10,12-octadecadien-1-yl)-D-glutaminy-L-threonyl-β-alanyl-D-tryptophyl-D-seryl-glycine, (6→2)-lactone  
**MF:** C<sub>46</sub>H<sub>66</sub>N<sub>8</sub>O<sub>11</sub>  
**FW:** 907.1  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/Acremonium sp.



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

### Laboratory Procedures

Aselacin C is supplied as a solid. A stock solution may be made by dissolving the aselacin C in the solvent of choice, which should be purged with an inert gas. Aselacin C is soluble in methanol and DMSO.

### Description

Aselacin C is a cyclic pentapeptolide antagonist of endothelin receptor type A (ET<sub>A</sub>) and ET<sub>B</sub> originally isolated from *Acremonium*.<sup>1</sup> It inhibits the binding of endothelin-1 to isolated bovine atrial membranes and porcine cerebral membranes, which endogenously express high levels of ET<sub>A</sub> and ET<sub>B</sub>, respectively, with IC<sub>50</sub> values of 60 and 80 µg/ml, respectively.

### Reference

1. Jackson, M., Burres, N.S., Karwowski, J.P., et al. Aselacins, novel compounds that inhibit binding of endothelin to its receptor. I. The producing organism, fermentation and biological activity. *J. Antibiot.* 47(5), 523-527 (1994).

#### 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, 07/07/2025

#### 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](http://WWW.CAYMANCHEM.COM)