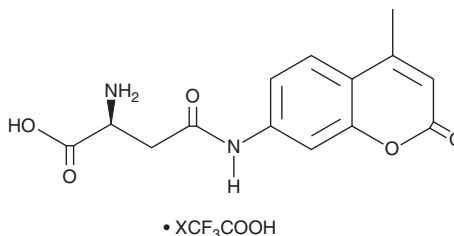


# PRODUCT INFORMATION

## L-Aspartic Acid $\beta$ -(7-amido-4-methylcoumarin) (trifluoroacetate salt) Item No. 45138

**Formal Name:** N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-asparagine, trifluoroacetate salt  
**MF:** C<sub>14</sub>H<sub>14</sub>N<sub>2</sub>O<sub>5</sub> • XCF<sub>3</sub>COOH  
**FW:** 290.3  
**Purity:**  $\geq$ 98%  
**Ex./Em. Max:** 360/440-450 nm  
**Supplied as:** A crystalline 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

L-Aspartic acid  $\beta$ -(7-amido-4-methylcoumarin) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the L-aspartic acid  $\beta$ -(7-amido-4-methylcoumarin) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. L-Aspartic acid  $\beta$ -(7-amido-4-methylcoumarin) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

L-Aspartic acid  $\beta$ -(7-amido-4-methylcoumarin) is a fluorescent probe for the detection of L-asparaginase activity.<sup>1</sup> It displays a weak fluorescent signal with an emission maximum of 400-405 nm upon excitation at 360 nm. Upon hydrolysis by L-asparaginase, 7-amido-4-methylcoumarin is released, displaying an increase in fluorescence intensity with an emission maximum of 440-450 nm upon excitation at 360 nm. It has been used to detect and quantify L-asparaginase activity in live cells and in serum.<sup>1,2</sup>

### References

1. Zlotnikov, I.D. and Kudryashova, E.V. A novel approach for the activity assessment of L-asparaginase formulations when dealing with complex biological samples. *Int. J. Mol. Sci.* **26(11)**, 5227 (2025).
2. Mononen, I.T., Kaartinen, V.M., and Williams, J.C. A fluorometric assay for glycosylasparaginase activity and detection of aspartylglycosaminuria. *Anal. Biochem.* **208(2)**, 372-374 (1993).

#### 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, 03/25/2026

#### 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