

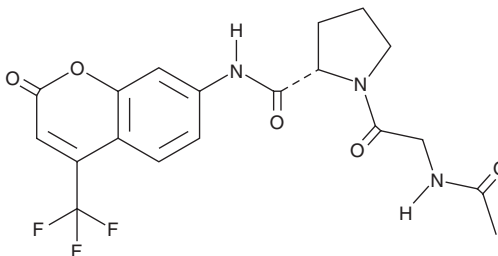
# PRODUCT INFORMATION



## Ac-Gly-Pro-AFC

Item No. 43373

**CAS Registry No.:** 886993-02-8  
**Formal Name:** N-acetylglycyl-N-[2-oxo-4-(trifluoromethyl)-2H-1-benzopyran-7-yl]-L-prolinamide  
**Synonym:** Ac-GP-AFC  
**MF:** C<sub>19</sub>H<sub>18</sub>F<sub>3</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 425.4  
**Purity:** ≥98%  
**Ex./Em.:** 400/505 nm  
**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

Ac-Gly-Pro-AFC is supplied as a solid. A stock solution may be made by dissolving the Ac-Gly-Pro-AFC in the solvent of choice, which should be purged with an inert gas. Ac-Gly-Pro-AFC is soluble (≥10 mg/ml) in ethanol and DMSO.

### Description

Ac-Gly-Pro-AFC is a fluorogenic substrate for fibroblast activation protein (FAP).<sup>1</sup> Upon enzymatic cleavage by FAP, 7-amino-4-trifluoromethylcoumarin (AFC) is released and its fluorescence can be used to quantify FAP activity. AFC displays excitation/emission maxima of 400/505 nm, respectively.

### Reference

1. Edosada, C.Y., Quan, C., Wiesmann, C., *et al.* Selective inhibition of fibroblast activation protein protease based on dipeptide substrate specificity. *J. Biol. Chem.* **281(11)**, 7437-7444 (2006).

#### 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, 06/26/2025

#### 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