

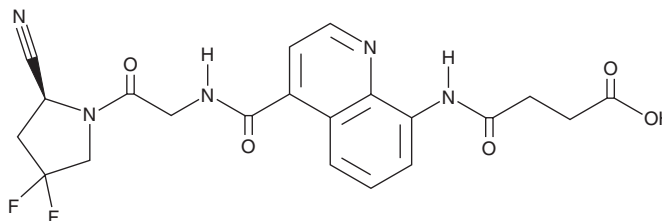
PRODUCT INFORMATION



OncoFAP

Item No. 44115

CAS Registry No.: 2639365-69-6
Formal Name: 4-[[4-[[[2-[(2S)-2-cyano-4,4-difluoro-1-pyrrolidinyl]-2-oxoethyl]amino]carbonyl]-8-quinoliny]amino]-4-oxo-butanoic acid
Synonym: Onco Fibroblast Activation Protein
MF: C₂₁H₁₉F₂N₅O₅
FW: 459.4
Purity: ≥95%
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

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

Description

OncoFAP is a fibroblast activation protein (FAP) ligand.¹ It binds to human FAP ($K_d = 0.68$ nM) and contains a carboxylic acid moiety that can be functionalized with various payloads for FAP-targeted delivery. OncoFAP has been conjugated to a variety of fluorophores and radioligands, which selectively accumulate in FAP-positive tumors in mice.

Reference

1. Millul, J., Bassi, G., Mock, J., *et al.* An ultra-high-affinity small organic ligand of fibroblast activation protein for tumor-targeting applications. *Proc. Natl. Acad. Sci. U S A* **118(16)**, e2101852118 (2021).

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, 01/20/2026

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