

# PRODUCT INFORMATION



## FOY-251

Item No. 36051

**CAS Registry No.:** 71079-09-9  
**Formal Name:** 4-[[4-[(aminoiminomethyl)amino]benzoyl]oxy]-benzeneacetic acid, monomethanesulfonate

**MF:** C<sub>16</sub>H<sub>15</sub>N<sub>3</sub>O<sub>4</sub> • CH<sub>3</sub>SO<sub>3</sub>H

**FW:** 409.4

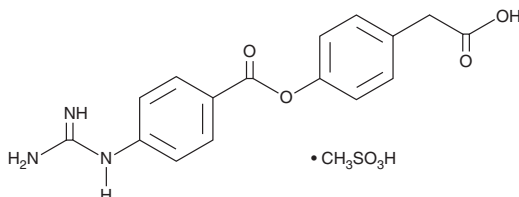
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 266 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

FOY-251 is supplied as a solid. A stock solution may be made by dissolving the FOY-251 in the solvent of choice, which should be purged with an inert gas. FOY-251 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of FOY-251 in these solvents is approximately 3 and 2 mg/ml, respectively.

### Description

FOY-251 is an active metabolite of the antiviral protease inhibitor camostat (Item No. 16018).<sup>1</sup> It inhibits transmembrane serine protease 2 (TMPRSS2; IC<sub>50</sub> = 33.3 nM).<sup>2</sup> FOY-251 inhibits entry of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) pseudotyped particles in Calu-3 cells (EC<sub>50</sub> = 178 nM) and reduces viral titers in SARS-CoV-2-infected Calu-3 cells when used at concentrations ranging from 1 to 10 μM.<sup>1</sup>

### References

1. Hoffmann, M., Hofmann-Winkler, H., Smith, J.C., *et al.* Camostat mesylate inhibits SARS-CoV-2 activation by TMPRSS2-related proteases and its metabolite GBPA exerts antiviral activity. *EBioMedicine* **65**, 103255 (2021).
2. Shrimp, J.H., Kales, S.C., Sanderson, P.E., *et al.* An enzymatic TMPRSS2 assay for assessment of clinical candidates and discovery of inhibitors as potential treatment of COVID-19. *bioRxiv* 167544 (2020).

#### 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, 12/19/2022

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