

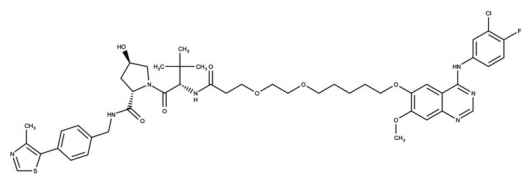
Gefitinib-based PROTAC 3

VHL-recruiting PROTAC targeting on EGFR

Gefitinib-based PROTAC 3 is targeting on EGFR, consisting of Gefitinib as the warhead linked to VHL-recruiting ligand.

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|---------------|---|
| Appearance | White solid. |
| CAS | 2230821-27-7 |
| Couple Target | EGFR, Ligase [E3] |
| Couple Type | Inhibitor, Ligand |
| Formula | $C_{47}H_{57}ClFN_7O_8S$ |
| MW | 934.5 |
| Purity | ≥99% |
| Solubility | Soluble in DMSO (75 mg/mL). Insoluble in water. |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)