

Product Information

Description	Gefitinib-based PROTAC 3 contains an EGFR binding element and a von Hippel-Lindau ligand joined by a linker which induces EGFR degradation.
Molecular Formula:	C ₄₇ H ₅₇ ClFN ₇ O ₈ S
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	934.5 Da
Cas No.	2230821-27-7

Chemical Structure

