

Product Information

Description	CFT7455 is an anti-cancer agent that binds with high affinity to the cereblon E3 ligase.
Molecular Formula:	$C_{28}H_{27}N_3O_4$
Supplied As:	Solid
Stability:	At least 6 months at $-20^{\circ}C$. Avoid freeze/thaw cycles.
Storage:	$-20^{\circ}C$
MW:	469.5 Da
Cas No.	2504235-67-8

Chemical Structure

