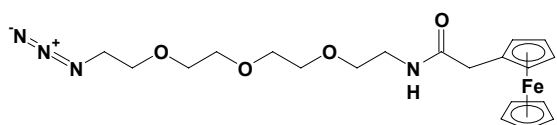


Ferrocene-PEG3- Azide

Cat. No.	Amount
CLK-CCA-9298-1	1 mg
CLK-CCA-9298-10	10 mg



Molecular Formula: C₂₀H₂₆FeN₄O₄

Molecular Weight: 442.29 g/mol

Storage conditions: store at -20 °C

Shelf life: 12 months

Appearance: yellow/brown, lyophilized powder/pellet

Purity: >90%(HPLC), +/- 2 daltons (ESI-MS)

For research use only!