

Na-Ascorbate

L-Ascorbic acid sodium salt

Cat. No.	Amount
CLK-MI005-10	10 mg
CLK-MI005-50	5 x 10 mg

CAS number: 134-03-2

Molecular formula: $C_6H_7O_6Na$

Molecular weight: 198.11 g/mol

Storage conditions: store at ambient temperature, store dry

Purity: > 99 %

Appearance: white powder

Melting point: 220 °C

Solubility: water

Shelf life: 12 months

For research use only!