

**Succinic Acid - 0.5 M**

any pH value
useful buffer range: pH 3.2 - 6.6
Butanedioic acid

Cat. No.	Amount
CSS-522	100 ml

Description:

Crystallization Stock Solution, 1M Succinic Acid adjusted with NaOH, sterile filtered

For in vitro use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: $\text{HOOCCH}_2\text{CH}_2\text{COOH}$

Molecular Weight: 118.09 g/mol

CAS#: 110-15-6

EC number: 203-740-4

Form: liquid

Concentration: 0.5 M

pH: arbitrary in 0.1 increments