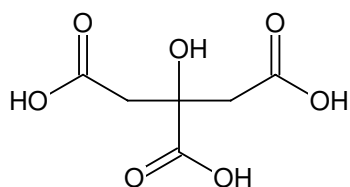


**Citric Acid - 1 M**

any pH value, useful buffer range: pH 2.2 - 6.5
sterile filtered

Cat. No.	Amount
CSS-508	100 ml



For *in vitro* use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: $\text{HOC}(\text{COOH})(\text{CH}_2\text{COOH})_2$

Molecular Weight: 192.12 g/mol

CAS#: 77-92-9

EC number: 201-069-1

Form: liquid

Concentration: 1 M

Description:

Crystallization stock solution: 1 M Citric Acid adjusted to any pH value within the buffer range pH 2.2 - 6.5 in 0.1 increments, adjusted with NaOH, also available without pH-adjustment

