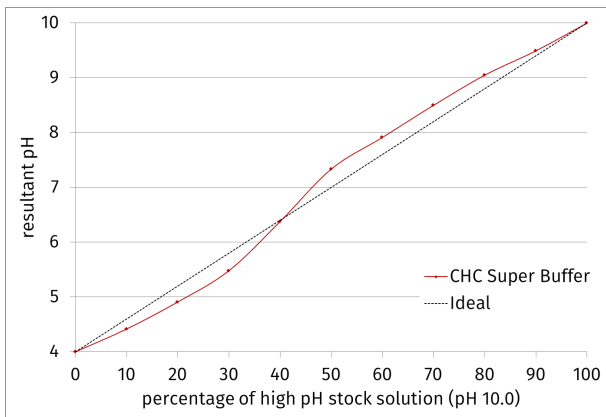




CHC pH 10.0 - 1 M

Buffer System: Citric Acid, HEPES, CHES
sterile filtered

Cat. No.	Amount
CSS-403	100 ml



CHC Super Buffer: Mixing low (CSS-402) and high (CSS-403) pH stock solutions results in an almost linear pH function

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula:

$C_6H_8O_7$
 $C_8H_{18}N_2O_4S$
 $C_8H_{17}NO_3S$

Molecular Weight: 192.12 g/mol; 238.31 g/mol; 207.29 g/mol

CAS#: 77-92-9, 7365-45-9, 103-47-9

Form: liquid

Concentration: 1 M

pH: 10.0, adjusted with NaOH

Description:

Crystallization Stock Solution - CHC Buffer System: Citric Acid, HEPES and CHES (2:3:4 Molar Ratio), pH 10.0 adjusted with NaOH

Citric Acid

Molecular Formula: $C_6H_8O_7$
Molecular Weight: 192.12 g/mol
CAS Number: 77-92-9

HEPES

Molecular Formula: $C_8H_{18}N_2O_4S$
Molecular Weight: 238.31 g/mol
CAS Number: 7365-45-9

CHES

Molecular Formula: $C_8H_{17}NO_3S$
Molecular Weight: 207.29 g/mol
CAS Number: 103-47-9



Related Products:

CHC pH 4.0, #CSS-402