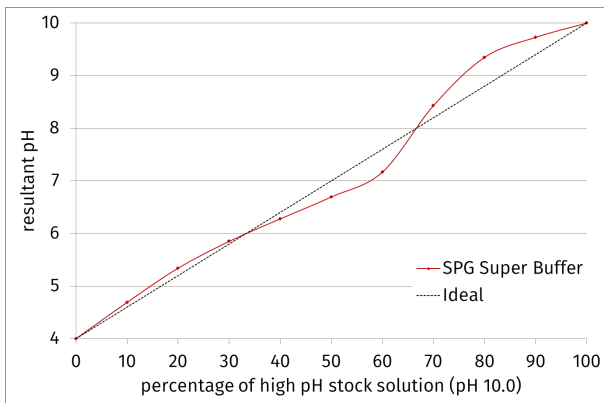




SPG pH 10.0 - 1 M

Buffer System: Succinic Acid, Sodium Dihydrogen Phosphate and Glycine
sterile filtered

Cat. No.	Amount
CSS-390	100 ml



SPG Super Buffer: Mixing low (CSS-389) and high (CSS-390) pH stock solutions results in an almost linear pH function

For *in vitro* use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula:

$C_4H_6O_4$

NaH_2PO_4

$C_2H_5NO_2$

Molecular Weight: 118.09 g/mol; 119.98 g/mol; 75.07 g/mol

CAS#: 110-15-6, 7558-80-7, 56-40-6

Form: liquid

Concentration: 1 M

pH: 10.0, adjusted with NaOH

Description:

Crystallization Stock Solution - SPG Buffer System: Succinic Acid, Sodium Dihydrogen Phosphate and Glycine (2:7:7 Molar Ratio), pH 10.0 adjusted with NaOH

Succinic Acid

Molecular Formula: $C_4H_6O_4$

Molecular Weight: 118.09

CAS Number: 110-15-6

Sodium Dihydrogen Phosphate

Molecular Formula: NaH_2PO_4

Molecular Weight: 119.98

CAS Number: 7558-80-7

Glycine

Molecular Formula: $C_2H_5NO_2$

Molecular Weight: 75.07

CAS Number: 56-40-6

Related Products:

SPG pH 4.0, #CSS-389