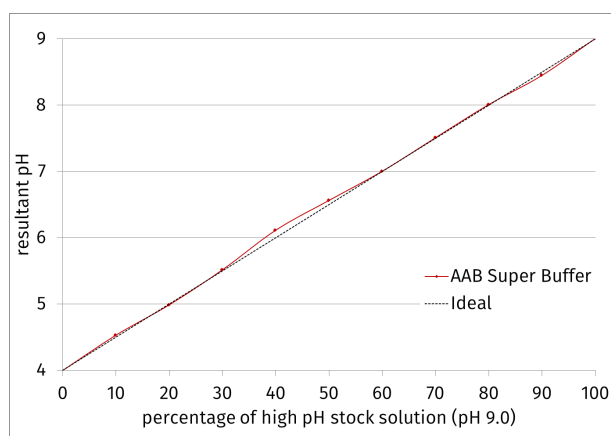


**AAB pH 9.0 - 1 M**

Buffer System: Sodium Acetate, ADA, Bicine
sterile filtered

Cat. No.	Amount
CSS-405	100 ml



AAB Super Buffer: Mixing low (CSS-404) and high (CSS-405) 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: C₂H₃NaO₂

C₆H₁₀N₂O₅

C₆H₁₃NO₄

Molecular Weight: 82.03 g/mol; 190.15 g/mol; 163.17 g/mol

CAS#: 127-09-3, 26239-55-4, 150-25-4

Form: liquid

Concentration: 1 M

pH: 9.0, adjusted with NaOH

Description:

Crystallization Stock Solution - AAB Buffer System: Sodium Acetate, ADA and Bicine (1:1:1 Molar Ratio), pH 9.0 adjusted with NaOH

Sodium Acetate

Molecular Formula: C₂H₃NaO₂

Molecular Weight: 82.03 g/mol

CAS Number: 127-09-3

ADA

Molecular Formula: C₆H₁₀N₂O₅

Molecular Weight: 190.15 g/mol

CAS Number: 26239-55-4

Bicine

Molecular Formula: C₆H₁₃NO₄

Molecular Weight: 163.17 g/mol

CAS Number: 150-25-4

Related Products:

AAB pH 4.0, #CSS-404