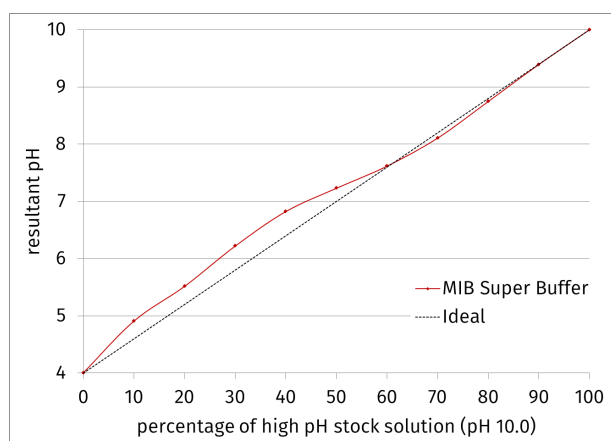


**MIB pH 10.0 - 1 M**

Buffer System: Malonic Acid, Imidazole, Boric Acid
sterile filtered

Cat. No.	Amount
CSS-401	4 x 25 ml



MIB Super Buffer: Mixing low (CSS-400) and high (CSS-401) 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:

$C_3H_4O_4$

$C_3H_4N_2$

H_3BO_3

Molecular Weight: 104.06 g/mol; 68.08 g/mol; 61.83 g/mol

CAS#: 141-82-2, 288-32-4, 10043-35-3

Form: liquid

Concentration: 1 M

pH: 10.0, adjusted with NaOH

Description:

Crystallization Stock Solution - MIB Buffer System: Malonic Acid, Imidazole and Boric Acid (2:3:3 Molar Ratio), pH 10.0 adjusted with NaOH

Malonic Acid

Molecular Formula: $C_3H_4O_4$

Molecular Weight: 104.06 g/mol

CAS Number: 141-82-2

Imidazole

Molecular Formula: $C_3H_4N_2$

Molecular Weight: 68.08 g/mol

CAS Number: 288-32-4

Boric Acid

Molecular Formula: H_3BO_3

Molecular Weight: 61.83 g/mol

CAS Number: 10043-35-3

**Related Products:**

MIB pH 4.0, #CSS-400