



Revised: June 24, 2024

Product Information

1X TAE (1L) Buffer Powder Packets (box of 50 packets)

Catalog Number: 22031

Unit Size: 50 packets (each packet makes 1 L of 1X buffer)

Storage and Handling

Store at room temperature. Product is stable for at least 12 months from date of receipt when stored as recommended. Handle with universal laboratory safety precautions.

Product Description

TAE Buffer (Tris-acetate-EDTA) is widely used for electrophoresis of nucleic acids in agarose and polyacrylamide gels. The buffer may be used as running buffer for gel electrophoresis, or when casting agarose gels. 1X TAE (1L) Buffer Powder Packets may be used as a running buffer for Precast GelRed® Agarose Gels (see Related Products).

Procedure

Prepare 1X TAE buffer by dissolving the contents of the buffer packet in 1 L of dH $_2$ O. The final composition of the 1X buffer will be 40 mM Tris-acetate, 1 mM EDTA, pH \sim 8 (do not adjust pH). 1X TAE buffer may be stored at room temperature indefinitely as long as the solution remains clear.

Note: We recommend using the same buffer for both casting and running gels for electrophoresis.

Related Products

Cat. No.	Product
41041	Precast GelRed® Agarose Gels, 1% Agarose/TAE
41003	GelRed® Nucleic Acid Gel Stain, 10,000X in Water
41001	GelRed® Nucleic Acid Gel Stain, 3X in Water
41011	GelRed® Prestain Plus 6X DNA Loading Dye
41029	GelRed® Agarose LE
41005	GelGreen® Nucleic Acid Gel Stain, 10,000X in Water
41030	GelGreen® Agarose LE
41032	EMBER500™ RNA Prestain Loading Dye
E90005	Gel-Bright™ Laser Diode Gel Illuminator
31030	DNA Gel Extraction Kit
31022	Ready-to-Use 1 kb DNA Ladder
31032	Ready-to-Use 100 bp DNA Ladder
99962-1	6X DNA Loading Buffer (Blue)
99859-1	6X DNA Loading Buffer (Orange)
31041	Forget-Me-Not™ EvaGreen® qPCR Master Mix (2-Color Tracking)
31000-T	EvaGreen® Dye, 20X in water (trial size), 1 mL
31077-T	EvaGreen® Plus Dye, 20X in water (trial size), 1 mL
31028	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit
41020	DNAzure® Blue Nucleic Acid Gel Stain
41028	Agarose LE, Ultrapure Molecular Biology Grade
41006	TBE Buffer, 5X (4L Cubitainer®)
41039	Go-Go™ Fast DNA Gel Running Buffer, 50X
22007	Activated Charcoal Decontamination Bags
41024-4L	Water, Ultrapure Molecular Biology Grade (4L Cubitainer®)

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Cubitainer is a registered trademark of Hedwin Corporation.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.