

Product Information

AccuClear® dsDNA Standard, 25 ng/uL

Catalog Number: 31029C

Unit Size: 1 mL

Storage and Handling

Store at 4°C. Product is stable for at least 12 months from date of receipt when stored as recommended. Use sterile, DNase-free pipettes, plates and tubes when handling the DNA standard to prevent degradation.

Product Description

The AccuClear® dsDNA Standard is a 25 ng/uL dsDNA standard, also provided in the AccuClear® Ultra High Sensitivity dsDNA Quantitation Kits. It contains dsDNA from calf thymus, 25 ng/uL in 10 mM Tris pH 7.5, 1 mM EDTA, 2 mM sodium azide. Potential applications include making custom DNA standard curves for quantitation.

Related Products

Catalog number	Product
31065	RNase-Free Calf Thymus DNA, 1 mg/mL
31006C	AccuBlue® High Sensitivity dsDNA Standards, Set of 8
31007C	AccuBlue® Broad Range dsDNA Standards, Set of 9
31006	AccuBlue® High Sensitivity dsDNA Quantitation Kit with DNA Standards
31008	AccuBlue® High Sensitivity dsDNA Quantitation Solution
31007	AccuBlue® Broad Range dsDNA Quantitation Kit with DNA Standards
31009	AccuBlue® Broad Range dsDNA Quantitation Solution
31028	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit with DNA Standards
31027	AccuClear® Ultra High Sensitivity dsDNA Quantitation Solution
31066	AccuGreen™ High Sensitivity dsDNA Quantitation Kit for Qubit®
31060	AccuBlue® NextGen dsDNA Quantitation Kit
41003	GelRed® Nucleic Acid Gel Stain, 10,000X in water
31000	EvaGreen® dye, 20X in water
31041	Forget-Me-Not™ EvaGreen® qPCR Master Mix
31043	Forget-Me-Not™ Universal Probe Master Mix
41006	TBE, 5X (4L)
41024	Water, Ultrapure Molecular Grade (4L)
22020	10X Phosphate Buffered Saline (PBS) (4L)
31022	Ready-to-Use 1 kb DNA Ladder
31032	Ready-to-Use 100 bp DNA ladder
31039	1 kb DNA Ladder in TE Buffer
31040	100 bp DNA Ladder in TE Buffer

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.

Qubit is a registered trademark of Thermo Fisher Scientific.