

ReadiUse™ PCR Reaction Buffer

Catalog Number: 17202

Unit Size: 50 mL

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

PCR (polymerase chain reaction) is a laboratory technique used to amplify DNA (deoxyribonucleic acid) segments. Our ready to use PCR buffer contains several components essential for the PCR reaction. The PCR reaction buffer helps to maintain a stable pH during the PCR reaction. It also contains a salt that helps to maintain the integrity of the DNA during the reaction and also enhances the activity of the DNA polymerase enzyme. MgCl₂ (magnesium chloride), a cofactor, is used for maintaining the activity of the DNA polymerase enzyme.