



NOTICE: ProZyme was purchased by Agilent in July 2018. Documents for products and product lots manufactured before August 2019 will contain references to ProZyme. For more information about these products and support, go to: www.agilent.com/en/contact-us.



PRODUCT DESCRIPTION

PRODUCT NAME: HORSERADISH PEROXIDASE CONJUGATE STABILIZER

PRODUCT CODE: CJ95

DESCRIPTION: Aqueous, protein-containing mixture preserved with 0.02% methylisothiazolone, 0.02% bromonitrodioxane and 20 ppm Proclin™ (Rohm and Haas)

STORAGE: Store at 4°C. Protect from direct exposure to light.

CJ95 HRP Conjugate Stabilizer is specifically designed for dilution of ProZyme Horseradish Peroxidase Conjugates. CJ95 HRP Conjugate Stabilizer prevents disassociation of the heme portion of the peroxidase molecule and maintains the conformation of the protein or antibody portion of the conjugate. It will not improve the stability of the bond between HRP and the antibody, protein or antigen.

DIRECTIONS FOR USE:

1. Use CJ95 Stabilizer directly from the bottle or dilute 1:1 in de-ionized water.
2. Dilute your conjugate using CJ95 Stabilizer to the optimum concentration for direct, sandwich or competitive assays.
3. Diluted conjugate is ready to use in your assay.
4. Store the diluted conjugate at 4°C. Protect from direct exposure to light.

NOTE: Phosphate buffers may cause precipitation at >10% total volume.