



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: ALKALINE PHOSPHATASE CONJUGATE STABILIZER

PRODUCT CODE: CJ90

DESCRIPTION: Aqueous, protein-containing mixture preserved with 0.02% methylisothiazolone and 0.02% bromonitrodioxane

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

CJ90 AP Conjugate Stabilizer is specifically designed for dilution of ProZyme Alkaline Phosphatase Conjugates. CJ90 AP Conjugate Stabilizer contains proprietary ingredients that prevent loss of catalytic activity and maintain conformation of the antibody, protein or antigen portion of the conjugate. It will not improve the stability of the bond between AP and the antibody, protein or antigen.

DIRECTIONS FOR USE:

1. Use CJ90 Stabilizer directly from the bottle or dilute 1:1 in de-ionized water.
2. Dilute your conjugate using CJ90 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.