

Product Information

Assay Buffer 1 (20X conc., 1.25mL, w/ DTT)

Product Number: **ASBF01**

Product Description

20X stock solution of assay buffer formulated for use in kinase assay applications.

Composition

400 mM HEPES, pH7.5 and 0.2 % TritonTM X-100, w/ 200 mM DTT

Prior to use

Dilute the buffer 20 times using ultrapure water and add DTT to achieve a final concentration of 2 mM. DTT is supplied as a 200 mM stock solution in a separate vial.

Storage and Stability

Store at -20°C.

Avoiding multiple freeze-thaw cycles is recommended to preserve product stability.