



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

CD45 (Human, recombinant, N-terminal His tag)
Catalog #: 30044

Description: Protein tyrosine phosphatase. Catalytic domain, (584-1256) (Gene bank accession number NM_080922) with N-terminal His tag, MW=75 kDa, expressed in a Baculovirus Sf9 expression system.

Source: Sf9 cells

Formulation: Recombinant proteins were eluted in 25 mM Tris-HCL, 75 mM NaCl, pH 8.0, 0.05% Tween, 5 mM DTT and 50% glycerol. Liquid at -20°C.

Stability: >12 months at -80°C

Purity: > 90% by SDS-PAGE

Specific Activity: 26 U/ug. One unit will hydrolyze 1 nmol p-nitrophenyl phosphate per minute at pH 7.4 and 30°C. Assay buffer: 50 mM HEPES, pH 7.4, 2 mM EDTA, 3 mM DTT, 100 mM NaCl, 50 mM pNPP. The specific activity of CD45 was determined using pNPP. Enzyme reaction condition: 20 mM pNPP, 2 min incubation at 30°C, 2 ug/ml enzyme.

Application: Useful for the study of enzyme kinetics, regulation, and to dephosphorylate target substrates.

References:

1. Rachmilewitz, J. et al., J. Biol. Chem. 278, 14059-14065
2. Fernandis, A.Z. et al., J. Biol. Chem. 278, 9536-9543

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com