

Data Sheet

Galactokinase

Human, recombinant, N-terminal His tag
Catalog #: 50275
 Lot# 121102-E2 Conc. 0.28 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 40 mM
 imidazole, 3 mM DTT, and 20% glycerol.

Stability: >12 months at -80 °C

References:

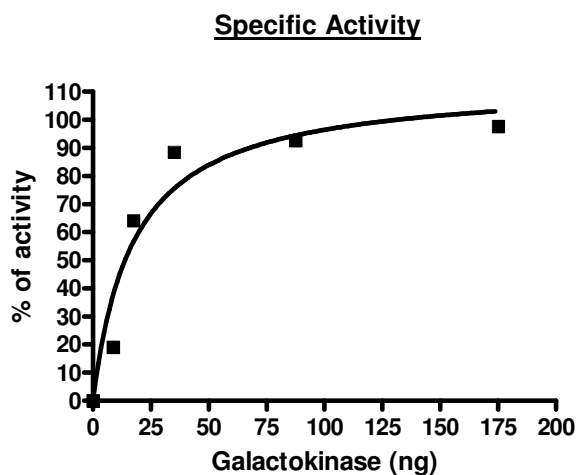
1. Thoden J.B. et al., *J. Biol. Chem.* **280** (10), 9662-9670 (2005).
2. Holden H.M. et al., *Cell. Mol. Life Sci.* **61** (19-20), 2471-2484 (2004).

Description: Human galactokinase (GenBank Accession No. NM_000154), full length with N-terminal His tag, MW=42 kDa, expressed in an *E. coli* expression system.

Specific activity: 102 pmole/min/μg
 A 50 μl Galactokinase reaction is conducted in a buffer containing 20 mM Tris (pH 7.4), 5 mM MgCl₂, 50 mM NaCl, 5 μM ATP, 1 mM DTT and 1 mM substrate D-Galactose at room temperature for 1 hour. ATP reduction is detected using Kinase-Glo[®] Luminescent Kinase Assay Platform.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2 μg Galactokinase

Lane 2:
 Protein Marker

MW: 42 kDa
Purity: ≥72%

