



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.



$\beta(1,4)$ -Galactosyltransferase

[$\beta(1,4)$ -Galactosyltransferase 1: B4GalT1]

SPECIFICATIONS

Product Code: GKT-GA14

Activity: ≥ 12 U/mg

Storage: -20°C

Shipped on ice pack for next day delivery.

Formulation: 50 mM Tris HCl
100 mM NaCl (pH 8.4)

$\beta(1,4)$ -Galactosyltransferase [$\beta(1,4)$ -Galactosyltransferase 1, EC 2.4.1.38] is a truncated version (amino acids 45-398) of human GalT, the enzyme responsible for the synthesis of Gal β 1-4GlcNAc. $\beta(1,4)$ -Galactosyltransferase was cloned and expressed in HEK 293F cells.

$\beta(1,4)$ -Galactosyltransferase (B4GalT1) transfers galactose from a donor substrate, UDP-galactose (UDP-Gal), to GlcNAc β 1-2Man units on glycoproteins and complex molecules.

Applications:

For *in vitro* galactosylation of glycoproteins such as monoclonal antibodies.

PRODUCT DESCRIPTION

Supplied Reagents:

WS0324 UDP-Gal
(3 X 10 mg; Uridine-5'-diphospho-galactose disodium salt)

WS0325 5x Reaction Buffer for GKT-GA14, 1 ml (50 mM MnCl₂, 500 mM MES, pH 6.5)

Molecular Weight: 39.5 kDa (by cDNA)

pH:

Recommended: 6.5
Range: 6.0 - 9.0

NOTE: A pH range of 6.0 to 9.0 may be used; at pHs below 6.5 enzyme activity may decrease; precipitation of MnCl₂ may occur at pHs above 7.0.

Stability: Store enzyme at -20°C . Avoid repeated freeze-thaw cycles as this decreases the efficacy of the enzyme. Dispense working aliquots (15 μg) after initial thaw if not utilizing entire quantity.

ASSAY

One unit of $\beta(1,4)$ -Galactosyltransferase is defined as the amount of enzyme required to release one mmole of UDP from UDP-Gal (measured as β NAD at 340 nm) per minute, at a dilution of ~0.1-0.2 U/ml, pH 8.0 and 30°C.

SUGGESTIONS FOR USE

Before use, briefly centrifuge the vial to ensure that all material is at the base of the vial. Ensure that reagents, substrates and laboratory-ware are free from contaminants and proteases.

The amount of enzyme required for galactosylation and the amount of galactosylation will vary depending on the target molecule (IgG, Fc-fusion, *etc.*) or application.

The recommended incubation time is 2-24 hours. Longer incubation times do not show intrinsic galactosidase activity. A time course to determine the optimal conditions for different intended targets is recommended.

The suggested buffer conditions for galactosylation are 10 mM MnCl₂, 100 mM MES, pH 6.5 with 10 mM UDP-Gal as the donor substrate.

Suggested Procedure for Galactosylation

To prepare a 500 μ g reaction using a standard IgG:

- 1) Prepare 1x Reaction Buffer.

For example, add 100 μ l of the supplied 5x Reaction Buffer to 400 μ l of ultrapure water.

The resulting 1x buffer contains 10 mM MnCl₂ with 100 mM MES, pH 6.5.

- 2) Prepare the target IgG to yield a final concentration of 10 mg/ml in 1x Reaction Buffer.
- 3) Dissolve one vial (10 mg) of UDP-Gal in 330 μ l of 1x Reaction Buffer.
- 4) Combine 500 μ g (50 μ l) of target IgG with 606 μ g (20 μ l) of UDP-Gal and 15 μ g (3 μ l) of $\beta(1,4)$ -Galactosyltransferase.
- 5) Adjust final volume to 100 μ l using 1x Reaction Buffer.
- 6) Incubate at 37°C for 7.5 hours; longer incubation times do not adversely affect enzyme activity.
- 7) Stop the incubation by freezing at -15 to -25°C.

NOTE: To avoid the potential for reaction-related artifacts affecting downstream analysis (e.g., in vitro and in vivo biological assays, PK/PD studies, mass spectrometric methods, etc.), the removal of excess reagents (e.g., by MWCO filtration, gel filtration, affinity chromatography, etc.) may be required.

