

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

Human UGT2B4 SUPERSOMES™

Catalog Number.....	456424	Storage Conditions..	STORE AT -80°C
Lot Number.....	2408296	Date Released	2024 September
		Expiration Date.....	2034 September

Package Contents.....0.5 mL
Protein Content.....5.0 mg/mL in 0.1M Tris pH 7.5
7-Hydroxy-4-trifluoromethylcoumarin
Glucuronidation Activity.....360 pmole/(min x mg protein)

PRODUCT DESCRIPTION: This activity is catalyzed by UGT2B4 which is expressed from human UGT2B4 cDNA using a baculovirus expression system. Baculovirus infected insect cells (BTI-TN-5B1-4) were used to prepare these membranes. A membrane preparation using wild type virus (Catalog No. 456400) should be used as a control for native activities.

ADVICE:

- Thaw rapidly in a 37°C water bath. Keep on ice until use.
- Aliquot to minimize freeze-thawing cycles. Less than 15% of the catalytic activity is lost after 8 freeze thaw cycles.
- Metabolite production using 7-hydroxy-4-trifluoromethylcoumarin glucuronidation as a substrate is linear with respect to enzyme concentration up to 0.8 mg/mL.
- Metabolite production with 7-hydroxy-4-trifluoromethylcoumarin is linear for at least 60 minutes (see graph). Other substrates may not exhibit similar linearity with respect to incubation time.

HAZARD WARNING: The product was produced using baculovirus (*Autographa californica*) infected insect cells (BTI-TN-5B1-4). This virus is not known to be pathogenic to humans or other mammals.

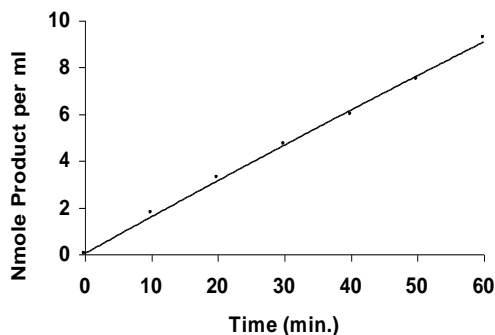
SAFETY INFORMATION: Safety assessment indicates this product is not hazardous, therefore no SDS (Safety Data Sheet) is provided. Use standard laboratory practices for the handling and disposal of Biosafety Level 1 materials.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

ASSAY METHOD: A 0.4 mL reaction mixture containing 0.25 mg/mL protein, 1 mM uridine diphosphoglucuronic acid (UDPGA), 10 mM magnesium chloride, 0.025 mg/mL alamethicin and 50 μ M 7-hydroxy-4-trifluoromethylcoumarin in 50 mM tris (pH 7.5) was incubated at 37°C for 20 minutes. After incubation, the reaction was stopped by the addition of 100 μ L 0.4 μ M labetalol in acetonitrile with 0.1% formic acid and centrifuged (10,000 x g) for 3 minutes. The product was detected by its Q1 and Q3 mass with negative polarity and quantitated by comparing the peak area ratio to a standard curve of 4-trifluoromethyl-7-hydroxycoumarin glucuronide.

Time Course of Product Formation



ANALYTICAL METHOD:

Materials

Column	2.1 x 50 mm 5 μ m C18 HPLC
Mobile Phase A	0.1% Formic Acid in dH ₂ O
Mobile Phase B	0.1% Formic Acid in Acetonitrile

Mass Transitions of MRM

Compound	Q1 Mass (amu)	Q3 Mass (amu)
Analyte- 4-trifluoromethyl-7-hydroxycoumarin glucuronide	405.1 \pm 0.2	288.7 \pm 0.2
Internal Standard- Labetalol	327.3 \pm 0.2	176.0 \pm 0.2

Gradient Separation Conditions

Time (minute)	Flow Composition of Mobile Phase A (%)	Flow Composition of Mobile Phase B (%)	Flow Rate (μ L /min)
0.0	95	5	750
2.5	5	95	750
2.6	95	5	750
3.6	95	5	750

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© 2013 DLS

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>



11 September 2024

Quality Assurance

Date