

β -Glucuronidase/Arylsulfatase

obtained from *Helix pomatia*

β -D-Glucuronoside glucuronosohydrolase, EC 3.2.1.31

Aryl-sulfate sulfohydrolase, EC 3.1.6.1

Cat. No. 10 127 060 001 2 ml

Cat. No. 10 127 698 001 10 ml

Version 09
Content version: March 2017

Store at +2 to +8°C

Product overview

Composition Enzyme mix of β -glucuronidase/arylsulfatase obtained from *Helix pomatia* in saline, stabilized with 0.02% sodium azide.

pH optimum

- The pH optimum value for β -glucuronidase activity is 4.5–5.0.
- The optimum for arylsulfatase activity is generally 6.2, but may be greater for some substrates in comparatively high concentrations; for instance, for 16.5 mM solutions of 2-hydroxy-5-nitrophenylhydrogen sulfate it is 7.2.

Inhibitors

- β -Glucuronidase activity is inhibited by:
 - D-glucuronic acid
 - D-galacturonic acid
 - D-glucaro-1,4-lactone (saccharolactone found in urine)
- Arylsulfatase activity is inhibited by phosphate.

Application

The enzyme preparation obtained from the Roman snail, *Helix pomatia*, which exhibits very strong β -glucuronidase and arylsulfatase activity, is widely used for the simultaneous hydrolysis of β -glucuronides (β -glucosiduronic acids) and sulfate esters in urine and other biological fluids (1, 2, 3, 4).

- Enzymatic hydrolysis of steroid β -glucuronides and sulfates
- Cell biology (removal of cell walls from yeasts in the preparation of protoplasts) (5)
- Enzyme immobilization studies (6)
- Determination of drugs in urine (7)

Storage/ Stability Undiluted, the aqueous solution of β -glucuronidase/arylsulfatase is stable at +2 to +8°C until the expiration date printed on the label.

Note: Aliquots portions of the diluted preparation may be stored at -15 to -25°C; they should not be thawed and refrozen more than once or twice; and storage at lower temperatures does not lengthen their expiration date beyond that of the product kept at +2 to +8°C.

Working concentration In many applications the product can be diluted with water immediately before use or used undiluted.

Note: This β -glucuronidase/arylsulfatase preparation is very concentrated and must be diluted for some applications. In particular for the preparation of protoplasts, the precise concentration to use for a given strain of yeast must be determined empirically.

Product description

Specificity of β -Glucuronidase The glycosides that β -D-glucuronic acid forms with a variety of compounds containing hydroxyl groups, hydrolyse readily in the presence of β -glucuronidase.

Such compounds include:

- steroids, such as estriol ($K_m = 0.42$ mM, pH 4.5), androsterone, pregnanediol, tetrahydrocortisone
- phenols, such as phenolphthalein ($K_m = 0.39$ mM), 4-nitro-phenol, 4-methylumbelliferone
- drugs, such as chloramphenicol and tetrahydrocannabinols
- metabolites, such as thyroxine and bilirubin

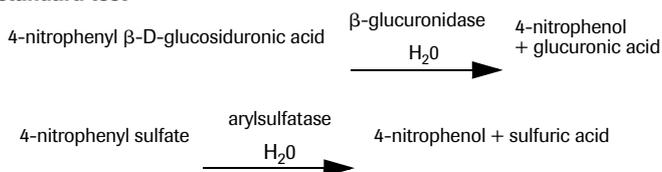
Polysaccharides that contain β -glucuronic acid residues, such as hyaluronic acid, are also hydrolyzed.

β -Glucuronidase is highly specific for the carbohydrate moiety: neither α -glucosides nor β -glucosiduronic acids are hydrolyzed.

Specificity of Arylsulfatase

Sulfate esters of many phenols are hydrolyzed in the presence of arylsulfatase. Examples are steroid sulfates, such as estronesulfate, 4-nitrophenyl hydrogen sulfate ($K_m = 1.8$ mM, pH 7.3), 4-nitro-pyrocatechol 2-sulfate ($K_m = 1.25$ mM, pH 7.5), and phenolphthalein disulfate.

Principle of the standard test



At a wavelength of 405 nm, the molar absorption coefficient of 4-nitrophenol is $1.85 \text{ mM}^{-1} \times 1 \times \text{cm}^{-1}$ at +25°C.

Steroids in urine

The various steroids found in urine may be present in one or more of three forms:

- (a) the free compound (in minor or trace quantities and amounts)
- (b) the sulfate (predominant in some cases)
- (c) the β -glucuronide (the predominant form in most cases)

Tab 1: Relative steroid proportions before hydrolysis are given in the following table:

Compound or category	free [%]	sulfate [%]	glucuronide [%]
7-Hydroxycorticosteroids	1	10–15	85–90
Pregnanediol	0	trace	≈ 100
Pregnanetriol	trace	trace	≈ 100
Estrone (O ₁)	1–3	10–15	85–89
Estradiols (O ₂)	1–3	5–10	90–95
Estriol (O ₃)	0–2	5–10	90–95
Androsterone	trace	20	80
Etiocholanolone	trace	10	90
Dehydroepiandrosterone (DHEA)	trace	≈ 100	trace
Epiandrosterone	trace	≈ 100	trace
11 β-Androsterone	trace	10	90
11 β-Etiocholanolone	trace	10	90
11 Ketoandrosterone	trace	trace	≈ 100
11 Ketoetiocholanolone	trace	trace	≈ 100

Methods for hydrolysis

Several methods of hydrolyzing steroid esters and glycosides are commonly used.

- For the sulfates of DHEA and androsterone, solvolysis is suitable; this involves treatment with excess organic solvent (*e.g.*, ethyl acetate, dioxan, or tetrahydrofuran) at a temperature of +38°C for 18–24 h.
- Acid hydrolysis at elevated temperatures is a more general method, but has two disadvantages: it can alter the structure and function of the steroids, and the resinified pigments formed need to be removed, because they are present in the extract.
- The third method, enzymatic hydrolysis (with β-glucuronidase and sulfatase), does not involve these drawbacks.

Procedure for the hydrolysis of glucuronides and sulfates in urine

Protocol

Please refer to the following table.

Step	Action
1	Adjust the pH of a portion of the sample (10 ml) to 5.5 by adding dilute acetic acid.
2	Add 1 ml acetate buffer (1 M pH 5.5) and 0.2 ml β-glucuronidase/ arylsulfatase solution.
3	Incubate at a temperature of +37°C for 16 h.
4	Cool and extract with an appropriate solvent (<i>e.g.</i> , chloroform or dichloromethane) to isolate the hydrolysis products.

Specific activity of Glucuronidase

Activity

Generally, the β-glucuronidase activity of the preparation is not as high with respect to steroid β-glucuronides, as values obtained from the hydrolysis of synthetic phenyl β-glucuronides indicate. However, under certain conditions, results obtained with phenolphthalein β-glucuronide may be comparable with those given by steroid glycosides, such as estradiol β-glucuronide. For instance, at +37°C and pH 4.5, a β-glucuronidase/arylsulfatase preparation that promotes the hydrolysis of 300 μmol of phenolphthalein β-glucuronide in 1 h, also promotes the hydrolysis of 441 μmol of estradiol β-glucuronide in 1 h.

Standard unit

The standard unit of β-glucuronidase activity is the enzyme activity that increases the rate of release of 4-nitrophenol from 4-nitrophenyl β-D-glucosiduronic acid at a temperature of +25°C and pH 4.5 by 1 μM.

Phenolphthalein unit

The phenolphthalein unit of β-glucuronidase activity is the enzyme activity that increases the rate of release of phenolphthalein from phenolphthalein β-D-glucosiduronic acid at a temperature of +38°C by 1 μM.

Approximately 4.5 standard units are equivalent to 5.5 phenolphthalein units.

Fishman unit

The Fishman unit (8) of β-glucuronidase activity is the enzyme activity that increases the rate of release of phenolphthalein from phenolphthalein β-D-glucosiduronic acid at a temperature of +38°C by 1 μg. Approximately 1 standard unit is equivalent to 22,000 Fishman units (1 phenolphthalein unit is equivalent to 19,000 Fishman units).

Specific β-glucuronidase activity

At +25°C and pH 4.5, the β-glucuronidase activity of 1 ml of the preparation is 4.5 standard units, equivalent to 5.5 phenolphthalein units or 100,000 Fishman units at +38°C.

Specific activity of Arylsulfatase

Standard unit

The standard unit of arylsulfatase activity is the enzyme activity that increases the rate of release of 4-nitrophenol from 4-nitrophenyl sulfate at a temperature of +25°C and pH 6.2 by 1 μM.

Phenolphthalein unit

The phenolphthalein unit of arylsulfatase activity is the enzyme activity that increases the rate of release of phenolphthalein from phenolphthalein disulfate at a temperature of +38°C and pH 6.2 by 1 μM.

Approximately 5.4 standard units are equivalent to 1 phenolphthalein unit.

Roy unit

The Roy unit of arylsulfatase activity is the enzyme activity that increases the rate of release of 4-nitroprocatechol from 2-hydroxy-5-nitrophenyl hydrogen sulfate (4-nitroprocatechol 2-sulfate) at a temperature of +38°C and pH 6.2 by 1 μg (9).

Approximately 1 standard unit is equivalent to 57,000 Roy units (1 phenolphthalein unit is equivalent to 308,000 Roy units).

Specific arylsulfatase activity

At +25°C and pH 6.2, the arylsulfatase activity of 1 ml of the preparation is 14 standard units, equivalent to 2.6 phenolphthalein units or 800,000 Roy units at +38°C.

Changes to previous version

Editorial changes

References

- 1 Bergmeyer, H.U., Grassel, M. & Walter, H.-E. (1983) in "Methods of Enzymatic Analysis" (Bergmeyer, H.U., eds.), 3. ed., Vol. 2, pp 206–207, VCH, Weinheim.
- 2 Stahl, P.D. & Fishman, W.H. (1984) in "Methods of Enzymatic Analysis", (Bergmeyer, H.U., eds.), 3. eds., Vol. 4, pp 246–256, VCH, Weinheim.
- 3 Tuite, M. et al. (1980) J. Biol. Chem. 255 8761.
- 4 Jarrige, P. Yon, J. & Hayle, M.F. (1963) Bull. Soc. Chim. Biol. 45 783.
- 5 von Hedenstroem, M. & Hoefer, M. (1974) Arch. Microbiol. 98 51.
- 6 Rapatz, E. et al. (1988) "Studies on the Immobilization of Glucuronidase (Part I and II)", in "Applied Biochemistry and Biotechnology", Vol. 19, pp 223–242.
- 7 Machata, G. & Kryspin-Eyner, K. (1970) Wiener klin. Wochenz. 47 849.
- 8 Wakabayashi, M. & Fishman, W.H. (1961) J. Biol. Chem 236 996.
- 9 Roy, A.B. (1970) in "Chemical and Biological Aspects of Steroid Conjugation" (Bernstein, S. & Solomon, S., eds.), pp 106, Springer, New York.

Trademarks

All other product names and trademarks are the property of their respective owners

Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

Disclaimer of License

For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

Contact and Support

To ask questions, solve problems, suggest enhancements and report new applications, please visit our [Online Technical Support Site](#).

To call, write, fax, or email us, visit sigma-aldrich.com, and select your home country. Country-specific contact information will be displayed.



Roche Diagnostics GmbH
Sandhofer Strasse 116
68305 Mannheim
Germany