

## User Guide

## Chorionic Gonadotropin Human

**C1063**

Storage Temperature -20 °C

CAS RN 9002-61-3

Synonyms: hCG, Choriogonin

## Product Description

pI: 2.95

Extinction Coefficient:<sup>1</sup>  $\epsilon^M = 1.41 \times 10^4$  (280 nm) $E^{1\%} = 3.6$  (280 nm)

Human chorionic gonadotropin (hCG) is a glycoprotein hormone produced by the chorionic tissue of the placenta. The glycoprotein hormone family includes luteinizing hormone (LH), follicle-stimulating hormone (FSH), and thyroid-stimulating hormone (TSH), in addition to hCG. The function of hCG is to maintain the corpus luteum and stimulate steroid secretion from the ovary in the beginning stages of gestation. hCG appears in the blood and urine during the first trimester of early pregnancy, and levels decrease thereafter. It has been used for superovulation in animals.<sup>2</sup>

The hCG consists of an  $\alpha$  subunit of 92 amino acids and a  $\beta$  subunit of 145 amino acids. The  $\alpha$  subunit is common among the family of glycoprotein hormones, whereas the hormone-specific  $\beta$  subunit, which exhibits different degrees of homology, may confer biologic specificity of the individual hormone. The amino acid sequences of the  $\alpha$  subunit<sup>4,5</sup> and the  $\beta$  subunit<sup>3,4</sup> and the crystal structure of Hcg<sup>6</sup> have been reported. NMR studies of the  $\alpha$  subunit have been reported.<sup>7,8</sup>

The molecular mass of hCG is ~37.9 kDa, with ~31% carbohydrate by weight. The theoretical molecular mass of 37.9 kDa is based on the native form, which contains 2 subunits. The  $\alpha$  subunit has a molecular mass of 14.9 kDa, with a polypeptide portion of ~10.2 kDa and a carbohydrate portion of ~4.7 kDa. The  $\alpha$  subunit is N-glycosylated at Asn<sup>52</sup> and Asn<sup>78</sup>.<sup>7</sup>

The  $\beta$  subunit has a molecular mass of 23 kDa, with a polypeptide portion of ~16.0 kDa and a carbohydrate portion of ~7.0 kDa.<sup>4,6,9,10</sup> The  $\beta$  subunit is N-glycosylated at Asn<sup>13</sup> and Asn<sup>30</sup>.<sup>11</sup> The structure and location of the O-glycosidic carbohydrate units attached to the  $\beta$  subunit of hCG, at Ser121, Ser127, Ser132, and Ser138, have been investigated.<sup>12</sup> The  $\beta$  subunit has been characterized by LC-MS.<sup>13</sup> A MALDI mass spectrometry report on the glycoforms of the hCG  $\beta$ -core fragment has been published.<sup>14</sup>

When hCG was used in combination with recombinant interferon- $\gamma$ , there was a significant cooperative induction of nitric oxide synthesis (iNOS) in a dose-dependent manner in mouse peritoneal macrophages. This suggests that hCG may provide a second signal for synergistic induction of NO synthesis.<sup>15</sup> The hCG has been used as a biomarker for such conditions as epithelial cancer<sup>16</sup> and germ cell tumors.<sup>17,18</sup>

The Chorionic Gonadotropin Human (C1063) is 0.2  $\mu$ m filtered. A vial contains ~2,500 IU. The calculated value is based on the USP method for human Chorionic Gonadotrophin to meet 2,500 IU per vial, with a tolerance of not less than 80.0% (2,000 IU) and not more than 125.0% (3,125 IU) of the potency stated.

## Preparation Instructions

When reconstituted with water to a concentration of 1,000 I.U./mL, the vial will also contain 0.04 M sodium phosphate buffer (pH ~7.2), and 40 mg/mL of mannitol. The hCG is also soluble in aqueous glycerol and glycols and is insoluble in ethanol.<sup>1</sup> Solutions should be sterile-filtered and not autoclaved.

## Precautions and Disclaimer

This product is for R&D use only, and not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

## Storage/Stability

Dilute aqueous solutions undergo rapid loss of activity when stored frozen, or heated, or if excess acid or base is added. Gelatin and serum proteins help to stabilize aqueous solutions of hCG. hCG is stable in a glycerol solution at 100 °C for one hour.<sup>19</sup>

Solutions reconstituted in water can be stored at -20 °C as single-use aliquots.

## References

1. Morgans, F.J. *et al.*, Endocrinology, 94(6), 1601-1606 (1974).
2. Hogan, B. *et al.*, Manipulating the Mouse Embryo, A Laboratory Manual, 2nd ed., Cold Spring Harbor Laboratory Press (Cold Spring Harbor, NY), p. 130 (1994).
3. Bellisario, R. *et al.*, J. Biol. Chem., 248(19), 6796-6809 (1973).
4. Morgan, F.J. *et al.*, J. Biol. Chem., 250(13), 5247- 5258 (1975).
5. Carlsen, R.B. *et al.*, J. Biol. Chem., 248(19), 6810- 6827 (1973).
6. Lapthorn, A.J. *et al.*, Nature, 369(6480), 455-461 (1994).
7. de Beer, T. *et al.*, FEBS Lett., 348, 1-6 (1994).
8. de Beer, T. *et al.*, Eur. J. Biochem., 241(1), 229- 242 (1996).
9. Bahl, O.P., J. Biol. Chem., 244(4), 567-574 (1969).
10. Swaminathan, N., and Bahl, O.P., Biochem. Biophys. Res. Commun., 40(2), 422-427 (1970).
11. Sisinni, L., and Landriscina, M., Adv. Exp. Med. Biol., 867, 159-176 (2015).
12. Kessler, M.J. *et al.*, J. Biol. Chem., 254(16), 7909- 7914 (1979).
13. Black, R.S., and Bowers, L.D., Methods Mol. Biol., 146, 337-354 (2000).
14. Jacoby, E.S. *et al.*, Clin. Chem., 46(11), 1796-1803 (2000).
15. Kim, H.M., and Moon, Y.H., Biochem. Biophys. Res. Commun., 229(2), 548-552 (1996).
16. Iles, R.K., Mol. Cell. Endocrinol., 260-262, 264-270 (2007).
17. Mitsios, J.V. *et al.*, Clin. Biochem., 47(7-8), 632-635 (2014).
18. Barlow, L.J. *et al.*, Nat. Rev. Urol., 7(11), 610-617 (2010).

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

## Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://www.sigmaaldrich.com/terms).

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://www.sigmaaldrich.com/offices).

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Merck and Sigma-Aldrich are trademarks of Merck or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2016-2025 Merck and/or its affiliates. All Rights Reserved.  
23251679 Rev 08/25

