

Product Information Sheet

Human CYP4F2BLR Bactosomes

Catalogue Number: CYP040

Product Description: Human CYP4F2 and human CYP-reductase coexpressed in Escherichia coli,

supplemented with purified human cytochrome b5

Amount: 1 nmol cytochrome P450

Storage Buffer: 50 mM Tris-acetate (pH 7.6), 250 mM sucrose, 0.25 mM EDTA

Storage: Store at -80°C. Avoid frequent temperature changes. Thaw on ice.

Intended Use: For laboratory (research) purposes only.

Batch Number:C4F2BLR001P450 Concentration:2.2 nmol/mlCytochrome b_5 Concentration:11.1 nmol/mlProtein Concentration:17.3 mg/ml

Specific P450 Content: 127 pmol/mg protein

Cytochrome *c* Reductase Activity: 208 nmol/min/mg protein

¹P450 Activity Data

Activity: Leukotriene B₄ ω-hydroxylase

²Linear with time up to: 30 min ³Linear with P450 up to: 60 pmol/ml

V_{max}: 1.7 pmol/min/pmol CYP

 K_m : 92 μ M

Approved by _____ Date ____14th December 2020

Notes

¹Unless otherwise stated, all assays are carried out at 37°C in 100 mM potassium phosphate buffer (pH 7.4) containing 5 mM magnesium chloride and an NADPH generating system [5 mM glucose 6-phosphate, 1 mM NADP⁺, 1 U/ml glucose 6-phosphate dehydrogenase].

²Determined at 20 µM substrate, using a P450 concentration of 50 pmol/ml.

Safety Data

The toxicological properties of this reagent have not been investigated. Exercise due care when handling.

IMPORTANT DISCLAIMER

- 1. Products supplied by Cypex may be harmful if misused. Any product ordered from Cypex must not be used for any purpose other than the intended use specified on the Product Information Sheet which accompanies the product when delivered. Please ensure that products supplied by Cypex are used safely, and, in particular, that the products do not come into direct human contact.
- 2. Cypex undertakes that all products supplied by it shall conform to the specification appearing on the Product Information Sheet which accompanies the product when delivered. In the event that products supplied by Cypex fail to meet such specification, Cypex will supply replacement products free of charge provided that the defective products have been returned to Cypex within a reasonable time of delivery (Cypex will reimburse reasonable costs incurred in returning defective products), and this shall be the only remedy available in respect of the supply of any defective products.
- 3. Cypex gives no undertakings or warranties in respect of products supplied other than the undertaking given at 2, above. Furthermore, Cypex has given and gives no undertaking or warranty nor does it make any representation, directly or indirectly, express or implied, about the quality, character or suitability of products supplied by Cypex.
- 4. Cypex shall have no liability in respect of loss or damage (whether direct or indirect or in the nature of loss of profit) suffered as a result of the supply of products by Cypex, or the subsequent use, storage or other application of the products, whether or not the products meet the specification referred to at 2, above.

³Determined at 20 µM substrate, using an incubation time of 10 min.