

POR Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # P242-30G

Lot # J425-4

Product Description

Recombinant full-length human POR was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC034277](#).

Gene Aliases

CYPOR

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

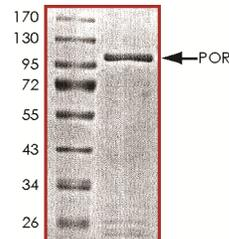
POR is a flavoprotein that donates electrons to all microsomal P450 enzymes, including the steroidogenic enzymes P450c17, P450c21, and CYP51A1. POR encodes an endoplasmic reticulum membrane oxidoreductase with an FAD-binding domain and a flavodoxin-like domain which binds two cofactors, FAD and FMN, which allow it to donate electrons directly from NADPH to all microsomal P450 enzymes (1). POR has been linked with various diseases, including apparent combined P450C17 and P450C21 deficiency, amenorrhea and disordered steroidogenesis, congenital adrenal hyperplasia and Antley-Bixler syndrome (2).

References

1. Shephard, E. A. et al: Isolation of a human cytochrome P-450 reductase cDNA clone and localization of the corresponding gene to chromosome 7q11.2. *Ann. Hum. Genet.* 53: 291-301, 1989.
2. Peterson, R. E. et al: Male pseudohermaphroditism due to multiple defects in steroid-biosynthetic microsomal mixed-function oxidases: a new variant of congenital adrenal hyperplasia. *New Eng. J. Med.* 313: 1182-1191, 1985.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity of POR was determined to be **>95%** by densitometry. Approx. MW **102 kDa**.

POR Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	P242-30G
Specific Lot Number	J425-4
Purity	>95%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.