

Purification of Recombinant Flavanone 3 β -Hydroxylase Using Bio-Scale CHT5-I Hydroxyapatite Column

Introduction

Flavanone 3 β -hydroxylase (FHT), a typical 2-oxoglutarate-dependent dioxygenase, catalyzes the 3 β -hydroxylation of 2S-flavanones to 2R,3R-dihydroflavonols which are intermediates in the biosynthesis of flavanoids, catechins, proanthocyanidins, and anthocyanidins in plants. To understand the complex reaction mechanism, the active site of the hydroxylase, and especially its interaction with the substrates and cofactors of the enzyme reaction, knowledge of the crystal structure of the protein is required.

The *E. coli*-expressed enzyme from *Petunia hybrida* was purified to apparent homogeneity by a three-step chromatographic procedure which included rapid, high resolution chromatography on a 5 ml Bio-Scale CHT5-I ceramic hydroxyapatite column (Bio-Rad, catalog # 751-0023).

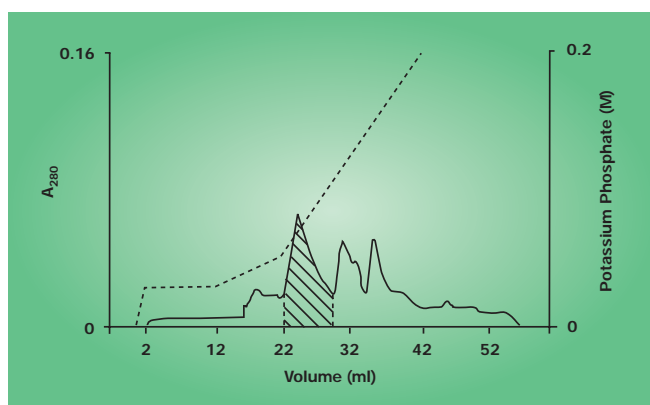


Fig. 1. Purification of recombinant flavanone 3 β -hydroxylase from *Petunia hybrida* on a Bio-Scale CHT5-I column. Protein load: 3 mg; Buffer A: 50 mM imidazole-HCl, 5 mM potassium phosphate, 2 mM DTT, pH 6.8; Buffer B: 200 mM potassium phosphate, 2 mM DTT, pH 6.8; Gradient: 12.5%–25% B in 10 ml, 25%–100% B in 40 ml; Flow rate: 1 ml/min; Detection: 280 nm.

Results

Recombinant flavanone 3 β -hydroxylase was expressed in *E. coli*. The cells were lysed and the resulting supernatant from a 45% saturated ammonium sulfate precipitation step was fractionated on a butyl HIC column. The FHT-containing fraction was concentrated and further fractionated using a gel filtration column. The FHT-containing fractions from the gel filtration column were pooled and loaded directly onto a 5 ml Bio-Scale CHT5-I column. Bound proteins were eluted by a two-step linear gradient from 25–50 mM in 10 ml and 50–200 mM potassium phosphate in 40 ml (Figure 1). Pure enzyme was obtained in the fractions containing 50–80 mM potassium phosphate. The purity of the enzyme was investigated by polyacrylamide gel electrophoresis (Figure 2).

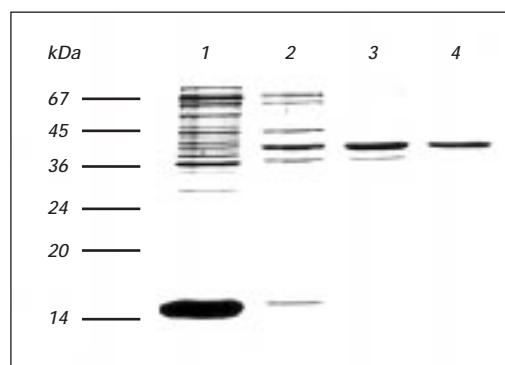
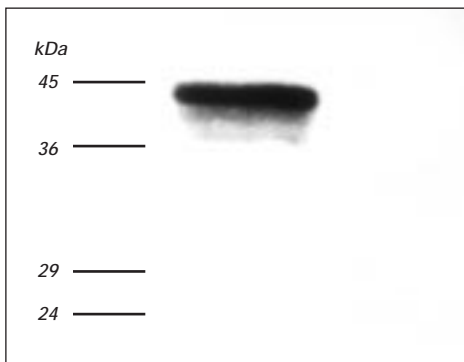


Fig. 2. SDS-PAGE analysis for *Petunia hybrida* flavanone 3 β -hydroxylase from *E. coli*. Fractions were run on a 12.5% Mini-PROTEAN[®] II gel and stained with Coomassie[®] blue stain. Lane 1: Soluble fraction of the *E. coli* cells expressing FHT (crude extract). Lane 2: Combined FHT-active fractions after hydrophobic interaction chromatography on Fractogel EMD Butyl (S). Lane 3: Combined FHT-active fractions after size exclusion chromatography on a Fractogel EMD BioSEC 650 (S). Lane 4: Purified FHT after chromatography on Bio-Scale CHT 5-I ceramic hydroxyapatite column.

Purification Table for Recombinant *Petunia hybrida* FHT from *E. coli* Cells (7 g)

PURIFICATION STEP	TOTAL PROTEIN (mg)	SPECIFIC ACTIVITY (U/mg)	PURIFICATION (FOLD)	TOTAL RECOVERY (%)
Crude Extract	236	0.05	1	100
Bio-Scale CHT5-I Column	0.535	2	40	9

Data is not provided for the butyl HIC and gel filtration steps



The purification procedure was monitored by immuno dot blot analysis for FHT protein (Figure 3) and enzyme activity assays of the individual fractions after each of the chromatographic steps. The table contains the essential purification parameters.

Data courtesy of Dr Richard Lukacin, Institute for Pharmaceutical Biology, University of Marburg, Germany
 Coomassie is a trademark of ICI.

Fig. 3. Immunoblot of FHT after Bio-Scale CHT5-I hydroxyapatite column probed with specific polyclonal antiserum from rabbit.



**Bio-Rad
Laboratories**

Life Science
Group

Website www.bio-rad.com U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771
 China 86-10-62051850/51 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France 01 43 90 46 90 Germany 089 318 84-0
 Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 Korea 82-2-3473-4460
 The Netherlands 0318-540666 New Zealand 64-9-4152280 Russia 7-95-4585822 Singapore 65-2729877 Spain (91) 661 70 85
 Sweden 46 (0)8 627 50 00 Switzerland 01-809 55 55 United Kingdom 0800-181134