

Purification of Functional Heterotrimeric G-Protein α -Subunits Using the Profinity eXact™ Fusion-Tag System on the Profinia™ Protein Purification System

Tony Ngo, Xiang Mao, and Kevin D. Ridge, Center for Membrane Biology, Department of Biochemistry and Molecular Biology, University of Texas Health Science Center, Houston, TX 77030

Introduction

Due to low natural abundance, many eukaryotic proteins require heterologous expression in order to obtain the milligram amounts of material often required for high-resolution structural studies. Prokaryotic expression systems have obvious advantages over eukaryotic cell expression systems, including better controlled growth in a variety of liquid media that can be supplemented with various labeled compounds to facilitate structure determination. A key limiting factor for prokaryotic expression systems is whether the eukaryotic protein of interest can be isolated in a pure and functional form. The use of affinity tags for isolation of heterologously expressed eukaryotic proteins has now become commonplace. Because the affinity tags can often interfere with the structure and/or function of the expressed protein, particularly in those cases where the amino- and/or carboxyl-terminus may contain a functionally important site(s), a cleavage site for a sequence specific protease (e.g. TEV protease, thrombin, etc.) is typically introduced between the protein of interest and the affinity tag. However, subsequent removal of the tag from milligram quantities of protein after affinity purification can be relatively costly, necessitating additional purification steps to remove the protease, and in many cases, may result in additional amino acids remaining on the protein of interest.

Here we describe the application of the Profinity eXact fusion-tag system (Oganesyan and Strong 2008; He et al. 2008), which is based on the interaction between the prodomain region of subtilisin BPN' and an active, mutant form of subtilisin BPN' (Ruan et al. 2004), integrated with the automated Profinia protein purification system to isolate milligram amounts of wild-type and mutant heterotrimeric G-protein α -subunits (G_{α}). The wild-type form of G_{α} , as well as a mutant (W254F), were found to be as active as G_{α} isolated from bovine retina.

Methods

Materials

Restriction endonucleases were purchased from New England Biolabs and Complete EDTA-free protease inhibitor cocktail tablets were obtained from Roche Applied Science. Phenylmethylsulfonyl fluoride (PMSF), isopropyl- β -D-thiogalactopyranoside (IPTG), and GDP were from Sigma-Aldrich, and bovine retinae were from W. Lawson Co. The Profinia purification system was from Bio-Rad Laboratories, Inc. A plasmid containing the gene for Chi6 (Skiba et al. 1996), an amino-terminal His6-tagged chimeric G_{α} protein that contains amino acids 1-215 and 295-350 from $G_{1\alpha}$ and an intervening sequence (amino acids 216-294) from $G_{11\alpha}$ was used as a source for the G_{α} protein gene used to make the expression construct. The pPAL5 expression vector, a fusion vector encoding the Profinity eXact tag (the modified 76 amino acid prodomain region of subtilisin BPN'), the 1 ml Bio-Scale™ Mini Profinity eXact cartridges containing immobilized S189 subtilisin BPN', and the 10 ml Bio-Scale Mini Bio-Gel® P-6 desalting columns were from Bio-Rad.

Insertion of the Chimeric G_{α} Gene Into pPAL5 and Mutant Construction

The construction of the Chi6 gene lacking the amino-terminal His6-tag and linker region has been described (Abdulaev et al. 2005a). The resulting gene was excised from the plasmid using HindIII and EcoRI and introduced into the pPAL5 vector using these same restriction sites. pPAL5 vector is essentially similar to the pPAL7 expression vector, a component of the Profinity eXact fusion-tag system, except for a few modifications in the multiple cloning sites downstream of the Profinity eXact fusion tag. Similarly, genes encoding the W207F and W254F mutants (Abdulaev et al. 2005a) were introduced into the pPAL5 vector using the HindIII and EcoRI restriction sites.

Expression of G_{α} Using the Profinity eXact Fusion-Tag System

E. coli BL21 (DE3) cells harboring a pPAL5/ G_{α} vector were grown in 200-250 ml of LB medium in the presence of 100 μ g/ml ampicillin at 26°C to $A_{550} \sim 0.3$, and then induced with 30 μ M IPTG for 12 hr at 26°C. The cell pellet was resuspended in 50 mM Tris-HCl, pH 8.0, containing 50 mM NaCl, 5 mM $MgCl_2$, 150 μ M GDP, 5 mM β -mercaptoethanol, 0.1 mM PMSF, and a protease inhibitor tablet and then disrupted by sonication.

Purification of G_{α} From the Profinity eXact Fusion-Tag Protein on an Automated Profinia System

The supernatant obtained by centrifugation of the cell lysate at 100,000 x g for 60 min was filtered through a 0.2 μ m filter, and then loaded onto a 1 ml cartridge packed with Profinity eXact purification resin integrated with a 10 ml Bio-Scale Mini Bio-Gel P-6 desalting cartridge on the Profinia protein purification system. All washing, elution, and column cleaning/regeneration steps were carried out using pre-programmed Bio-Rad methods and recommended buffers except that the binding and wash buffers also contained 5 mM $MgCl_2$, 150 μ M GDP, and 5 mM β -mercaptoethanol. The pre-elution incubation step with 100 mM NaF was for 2 hr. For the W207F and W254F mutant G_{α} constructs, purification using the Profinity eXact purification method was performed in tandem on the Profinia system, under the same conditions as described above for wild-type G_{α} .

Other Methods

The retinal heterotrimeric G-protein transducin (G_t) was isolated from bovine retina using a standard protocol (Fung et al. 1981), and $G_{t\alpha}$ separated from the $G_{t\beta\gamma}$ subunits as described (Abdulaev et al. 2005b). The intrinsic tryptophan fluorescence of the various G_{α} subunits in the absence and presence of aluminum fluoride (AlF_4^-) was measured essentially as described (Abdulaev et al. 2005b). Protein samples were analyzed by SDS-PAGE using Criterion™ XT Bis-Tris gels 4-12% (Bio-Rad) and XT MES running buffer (Bio-Rad) and visualized by Coomassie Blue staining. Protein determinations were done using the method of Peterson (Peterson 1977) with BSA as the standard.

Results and Discussion

Expression and Purification of Profinity eXact Fusion-Tagged G_{α} Subunits

Wild-type and two mutant forms (W207F and W254F) of a chimeric G_{α} subunit were expressed in *E. coli* as a Profinity eXact fusion-tag and purified on Profinity eXact purification resin integrated with a Bio-Scale Mini Bio-Gel P-6 desalting cartridge on the Profinia protein purification system. Figures 1A and 1B show a representative chromatogram and results from SDS-PAGE analysis, respectively, for wild-type G_{α} . The chromatogram shows that G_{α} can be effectively eluted from the Profinity eXact resin and desalted on the Bio-Scale Mini Bio-Gel P-6 cartridges with a good yield. Consistent with earlier G_{α} chimera expression studies (Abdulaev et al. 2005a), results using the Profinity eXact fusion-tag encoded by the

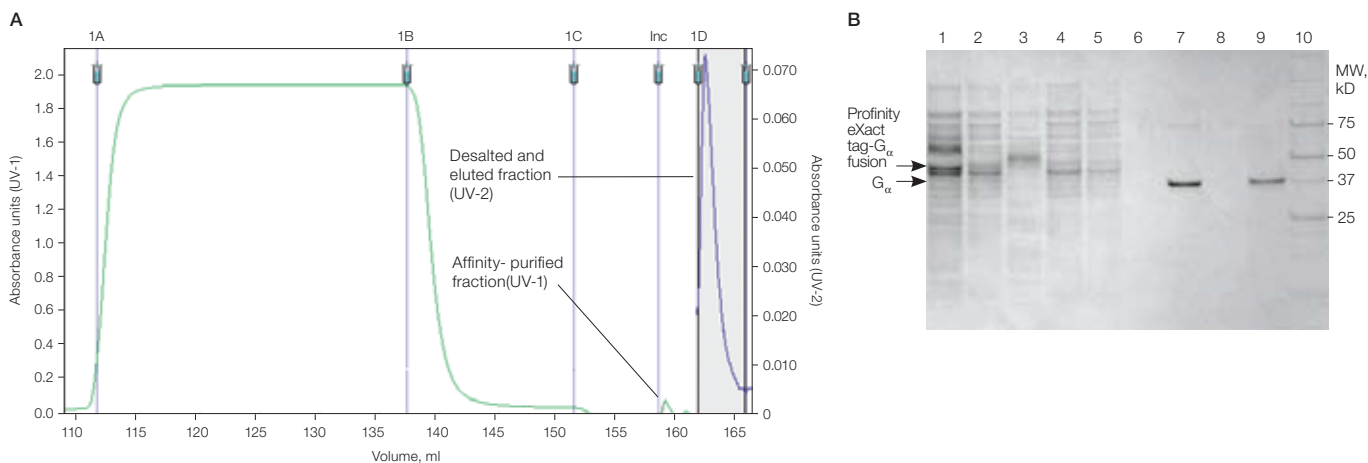


Fig. 1. G_{α} purification using Profinity eXact purification plus integrated desalting method on the Profinia system. **A**, chromatogram of Profinity eXact-tag- G_{α} fusion purification using Profinity eXact purification resin and a Bio-Scale Mini Bio-Gel P-6 desalting cartridge. The soluble fraction of the cell lysate was loaded onto the 1 ml Bio-Scale Mini Profinity eXact cartridge (1A, flowthrough) and the column was washed with 14 column volumes (CVs) of binding/wash 1 buffer, 100 mM sodium phosphate, pH 7.2 (1B, wash 1), followed by 6 CVs of wash 2 buffer, 100 mM sodium phosphate, 300 mM sodium acetate, pH 7.2 (1C, wash 2). Both buffers were supplemented with 5 mM $MgCl_2$, 150 μ M GDP, and 5 mM β -mercaptoethanol. The Bio-Scale Mini Profinity eXact cartridge was primed with 1 CV of elution buffer, 100 mM sodium phosphate, 100 mM sodium fluoride, pH 7.2, supplemented with 150 μ M GDP and 5 mM β -mercaptoethanol, and allowed to incubate for 2 hours (Inc, enzyme cleavage). G_{α} obtained by fluoride-triggered cleavage from the bound Profinity eXact tag, was eluted in 3.5 CVs of the same buffer and desalted using a 10 ml Bio-Scale Mini Bio-Gel P-6 cartridge (1D, elution). (—) peaks detected by UV1 monitor. (—) detected by UV2 monitor; **B**, SDS-PAGE analysis of fractions obtained at various steps of the purification process. Lane 1, total cell extract from IPTG-induced cells; 2, 3, soluble and insoluble extracts respectively, from IPTG-induced cells; 4, column flowthrough (1A); 5, column wash 1 (1B); 6, column wash 2 (1C); 7, elution fraction following desalting (1D); 8, blank; 9, G_{α} from bovine retina; and 10, Precision Plus Protein™ standards. Protein bands were visualized using Coomassie Blue stain and the positions of some of the molecular mass standards are shown on the right. The arrows on the left indicate the position of the Profinity eXact tag- G_{α} fusion and tag-free G_{α} .

pPAL5 vector clearly show that a majority of the fusion protein is present in the supernatant of the cell lysate (Figure 1B, compare lanes 1-3). Further, the soluble fraction containing the fusion protein binds to the Profinity eXact purification resin (Figure 1B, lane 4) and after two rounds of washing to remove contaminating proteins (Figure 1B, lanes 5 and 6) is subsequently cleaved from the Profinity eXact fusion-tag by the immobilized protease in the presence of 100 mM NaF and desalted on the integrated Bio-Scale Mini Bio-Gel P-6 desalting cartridge on the Profinia system. The resulting protein, which we refer to as ChiT (Abdulaev et al. 2005a), is present in a highly purified form (Figure 1B, lane 7) and co-migrates with $G_{t\alpha}$ isolated from bovine retina (Figure 1B, lane 9). Similar results were obtained for the W207F and W254F mutant forms of G_{α} when expressed and purified using the Profinity eXact fusion-tag system on the Profinia system (data not shown).

AlF_4^- -Dependent Tryptophan Fluorescence Changes of G_{α}

Adduct formation with AlF_4^- and the GDP/ Mg^{2+} bound form of G_{α} results in an increase in the intrinsic tryptophan fluorescence of the protein. In $G_{t\alpha}$, this increase in fluorescence has been largely attributed to a change in the environment of Trp-207 (Faurobert et al. 1993). As shown in Figure 2A, $G_{t\alpha}$ shows a ~30% increase in intrinsic fluorescence upon AlF_4^- treatment. Similarly, ChiT also shows ~30% increase in fluorescence emission upon AlF_4^- adduct formation (Figure 2B), suggesting that G_{α} prepared using the Profinity eXact fusion-tag system on the Profinia system is as functionally active as native $G_{t\alpha}$ isolated from bovine retina. Further, we have previously shown that G_{α} subunits expressed and purified using similar approaches are able to form stable heterotrimers with $G_{t\beta\gamma}$ subunits and undergo receptor-catalyzed guanine nucleotide exchange (Abdulaev et al. 2005b; Ridge et al. 2006a).

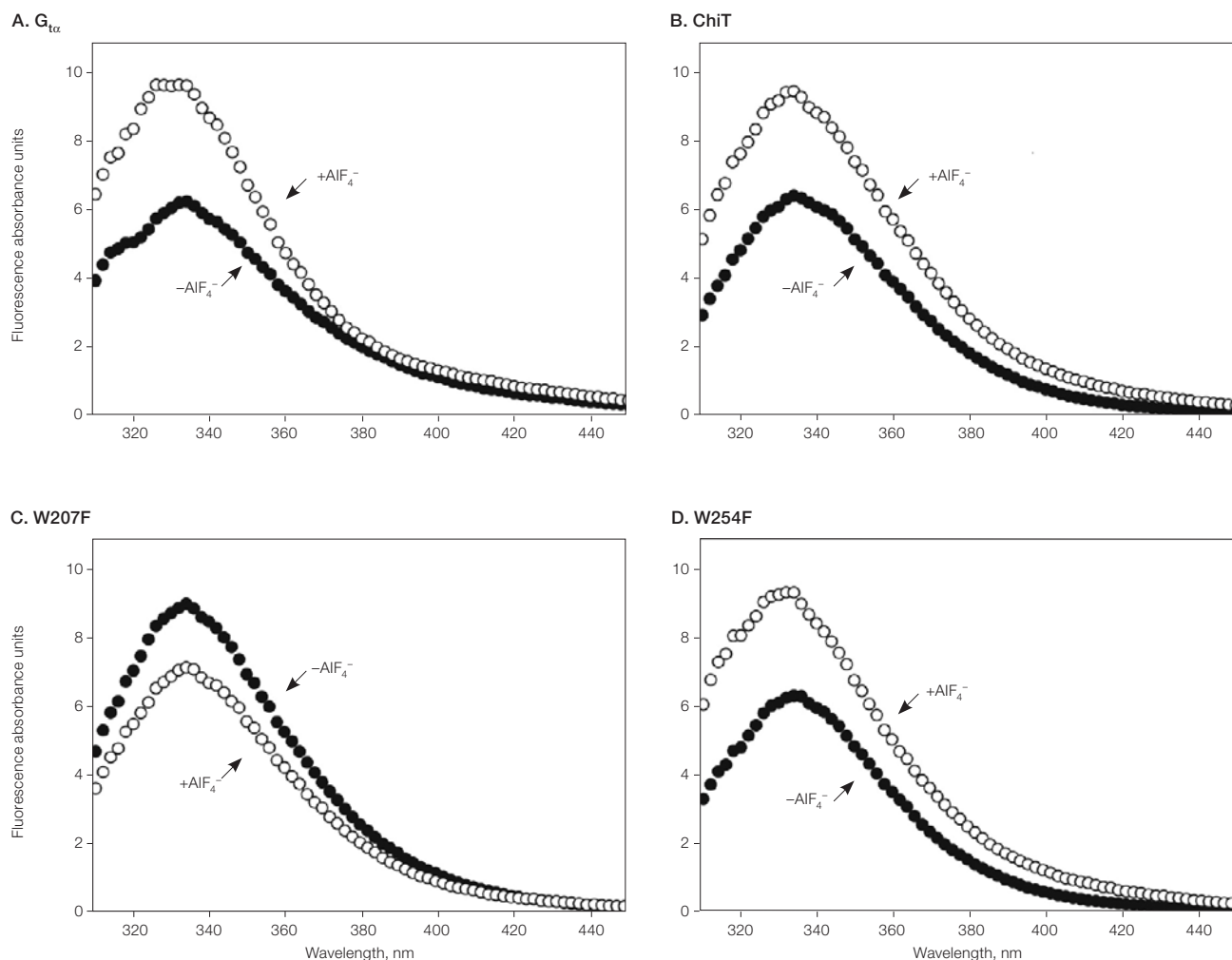


Fig. 2. Fluorescence changes accompanying AlF_4^- treatment of the GDP/ Mg^{2+} -bound form of wild-type and mutant forms of G_{α} isolated using the Profinity eXact fusion-tag purification method on the Profinia system. A, B, C, and D, represent emission spectra for $G_{t\alpha}$, ChiT, W207F G_{α} , and W254F G_{α} respectively. (●), fluorescence before and (○), fluorescence after the addition of AlF_4^- .

As expected, the W207F mutant did not show the corresponding fluorescence increase upon addition of AlF_4^- , but in fact showed a ~10% decrease (Fig. 2C). This may be attributed to a change in the environment of W254F. In this respect, the W254F mutant showed a similar increase in fluorescence upon addition of AlF_4^- as $\text{G}_{\alpha\text{T}}$ and ChiT.

Summary

We have shown using the Profinity eXact fusion-tag system integrated with the Profinia protein purification system that milligram quantities of a highly purified wild-type G protein α -subunit chimera (ChiT) as well as mutant forms of the protein can be isolated in a relatively quick and reproducible fashion. Further, purified ChiT and the W254F mutant of G_{α} are as functionally active as G_{α} isolated from natural sources. This technical advance should further facilitate efforts using high-resolution structural methods to investigate the conformation and dynamics of G_{α} in various states (Abdulaev et al. 2005a; Abdulaev et al. 2005b; Ridge et al. 2006a; Ridge et al. 2006b; Abdulaev et al. 2006).

Acknowledgment

Work in the Ridge laboratory on heterotrimeric G proteins is supported by NIH Grant EY016493.

References

- Abdulaev NG et al. (2005a). Bacterial expression and one-step purification of an isotope-labeled heterotrimeric G-protein α -subunit. *J Biol Chem* 280, 31-40.
- Abdulaev NG et al. (2005b). Heterotrimeric G-protein α -subunit adopts a "preactivated" conformation when associated with $\beta\gamma$ -subunits. *J Biol Chem* 280, 38071-38080.
- Abdulaev NG et al. (2006). The receptor-bound "empty pocket" state of the heterotrimeric G-protein α -subunit is conformationally dynamic. *Biochemistry* 45, 12986-12997.
- Faurobert E et al. (1993). Tryptophan W207 in transducin $\text{T}\alpha$ is the fluorescence sensor of the G protein activation switch and is involved in the effector binding. *EMBO J* 12, 4191-4198.
- Fung BK et al. (1981). Flow of information in the light-triggered cyclic nucleotide cascade of vision. *Proc Natl Acad Sci U S A* 78, 152-156.
- He X et al. (2008). Purification of tag-free recombinant proteins using the Profinity eXact fusion-tag system. *Bio-Rad Bulletin* 5656.
- Oganesyan N and Strong W (2008). Profinity eXact fusion-tag system performs on-column cleavage and yields pure native protein from lysate in less than an hour. *Bio-Rad Bulletin* 5652.
- Peterson GL (1977). A simplification of the protein assay method of Lowry et al. which is more generally applicable. *Anal Biochem* 83, 346-356.
- Ridge KD et al. (2006a). Conformational changes associated with receptor-stimulated guanine nucleotide exchange in a heterotrimeric G-protein α -subunit: NMR analysis of GTP γ S-bound states. *J Biol Chem* 281, 7635-7648.
- Ridge KD et al. (2006b). NMR analysis of rhodopsin-transducin interactions. *Vision Res* 46, 4482-4492.
- Ruan B et al. (2004). Engineering subtilisin into a fluoride-triggered processing protease useful for one-step protein purification. *Biochemistry* 43, 14539-14546.
- Skiba NP et al. (1996). Mapping of effector binding sites of transducin α -subunit using $\text{G}\alpha\text{T}/\text{G}\alpha\text{1}$ chimeras. *J Biol Chem* 271, 413-424.

Coomassie Blue is a trademark of BASF Aktiengesellschaft.

Complete is a trademark of Roche Applied Science.

Profinity eXact vectors, tags, and resins are exclusively licensed under patent rights of Potomac Affinity Proteins. This product is intended for research purposes only. For commercial applications or manufacturing using these products, commercial licenses can be obtained by contacting the Life Science Group Chromatography Marketing Manager, Bio-Rad Laboratories, Inc., 6000 Alfred Nobel Drive, Hercules, CA 94547, Tel (800)4BIORAD.

The composition and/or use of the T7 expression system is claimed in one or more patents licensed to Bio-Rad by Brookhaven Science Associates, LLC. A separate license is required for any commercial use, including use of these materials for research or production purposes by any commercial entity.

Information in this tech note was current as of the date of writing (2008) and not necessarily the date this version (rev A, 2008) was published.



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000