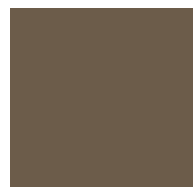




**Profinity eXact™**  
Fusion-Tag System



Affinity Tag Purification and Tag Removal



## Fast Forward to Tag-Free Protein

Protein fusion technology has greatly simplified the process of purifying recombinant proteins. Recombinant technology has made the expression of large numbers of proteins possible, usually in tagged format, allowing their use in standard chromatography applications.

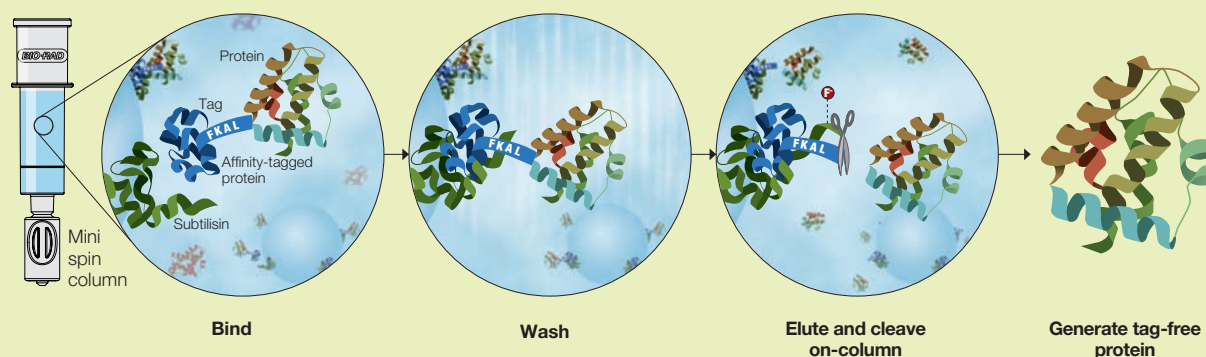
The Profinity eXact fusion-tag system addresses the bottlenecks in the tag-removal process during affinity-tagged protein purification. The Profinity eXact system is a novel *E. coli*-based system for the expression, detection, purification, and on-column cleavage of affinity-tagged proteins, without the addition of protease. The system offers single-step purification without the hassle and expense of cleavage enzymes, incubation times, or removal of reagents.

### Key Features

- Purification and processing of fusion-tagged proteins in a single step
- On-column cleavage in as little as 30 minutes
- Elimination of protease-addition step
- Precise cleavage at the N-terminus to generate native protein sequence



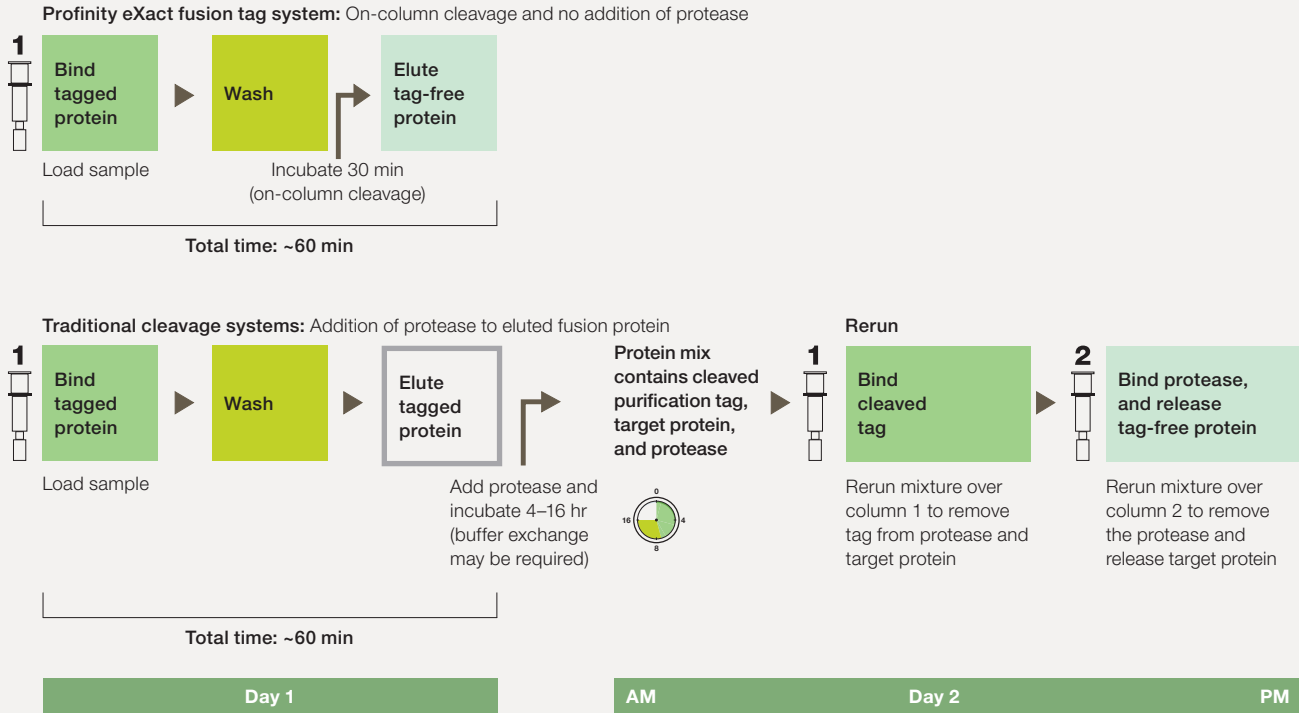
Profinity eXact fusion-tag system consists of expression vectors, chemi-competent cells, SOC growth media, spin columns, purification resin, cartridges, and detection reagents.



**Purification and on-column tag cleavage using the Profinity eXact fusion-tag system.** During sample application, immobilized subtilisin protease (ligand) recognizes and binds the affinity-tagged protein. Washing the column removes unbound contaminants. Application of fluoride-containing elution buffer triggers subtilisin to quickly and precisely cleave the tag from the fusion protein after the cleavage recognition sequence ending in FKAL. The tag remains tightly bound to the resin, and a highly purified protein with only its native amino acid sequence is released.

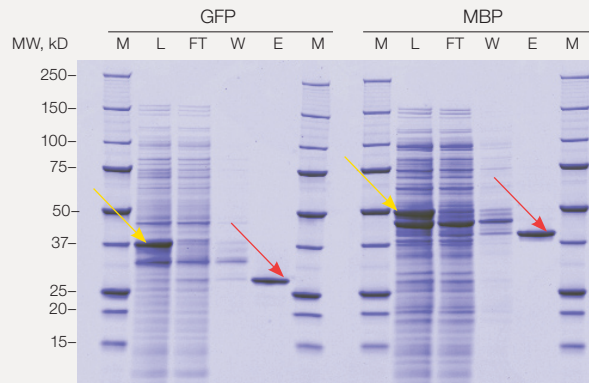
### One-Step Purification and Tag Cleavage

In experiments requiring tag cleavage, protease and tag removal problems are unavoidable. The Profinity eXact fusion-tag system easily overcomes these challenges with parallel purification and on-column cleavage. Availability of immobilized protease in the column appreciably shortens the purification process.



### Precise Cleavage and High Purity

Immobilized protease on a chromatography support is unique to Profinity eXact purification resin, distinguishing it from traditional methods of tag cleavage. Purity of the eluted protein is typically higher than for other affinity tag systems due to the specific recognition of the immobilized subtilisin protease for its tag sequence and the high fidelity of the protease reaction.



**SDS-PAGE analysis of proteins, GFP (26 kD) and MBP (40 kD), fused to Profinity eXact purification resin and purified by Bio-Scale™ Mini Profinity eXact cartridges.** Crude *E. coli* lysate (2 ml) containing the expressed proteins was loaded onto a 1 ml Bio-Scale Mini Profinity eXact cartridge on a BioLogic DuoFlow™ system with binding/wash buffer at 1 ml/min. The cartridge was washed with 10 column volumes (CV) of the same buffer at 1 ml/min. Proteins were eluted with 3 CV potassium fluoride buffer at 0.1 ml/min for 30 min at room temperature. Total purification time required to generate tag-free proteins without addition of protease was approximately 60 min. M, markers (Precision Plus Protein™ standards); L, load (crude lysate); FT, flowthrough; W, wash; E, eluate. (→), tagged proteins; (→), tag-free purified proteins.

