

Recombinant 6H-PFO* (Y181A,C459A)

Project#: Codex10121719	Lot#: 112217G	Concentration: 2.0mg/ml
--------------------------------	----------------------	--------------------------------

Description The coding region of PFO* (Y181A,C459A) was subcloned into pET28a vector with a 6-His tag at the N-terminus. Recombinant 6H-PFO* was expressed in BL21(DE3) host cells, induced by IPTG and purified from soluble lysate by Ni-sepharose affinity column.

MRGSHHHHHHGMASMTGGQOMGRDLYDDDDKDPSSRSKDIITDKNQSIDSG
ISSLSYNRNEVLASNGDKIESFVPKEGKKTGNKFIVVERQKRSLTTSPVD
ISIIDSVNDRITYPGALQLADKAFVENRPTILMVKRKPININIDLPLGLKGE
NSIKVDDPTYGKVS GAIDELVSKWNEKYSSTHTLPARTQASESMVYSKSQ
ISSALNVNAKVLENSLGVDFAVANNEKVMILAYKQIFYTVSADLPKNP
SDLFDDSVTFNDLKQKGVSNAPPLMVSNAVAYGRTIYVKLETTSSSKDVQ
AAFKALIKNTDIKNSQQYKDIYENSSTAVVLGGDAQEHNKVVTKDFDEI
RKVIKDNATFSTKNPAYPISYTSVFLKDNSVAAVHNKTDYIETTSTEYSK
GKINLDHSGAYVAQFEVAWDEVSYDKEGNEVLTHKTTWDGNYQDKTAHYST
VIPLEANARNIRIKAREATGLAWEWWRDVISEYDVPLTNNINVS IWGTTL
YPGSSITYN

Source *E. coli.*

Molecular Mass Approximately 60kDa.

Purity ≥90%, as determined by SDS-PAGE.

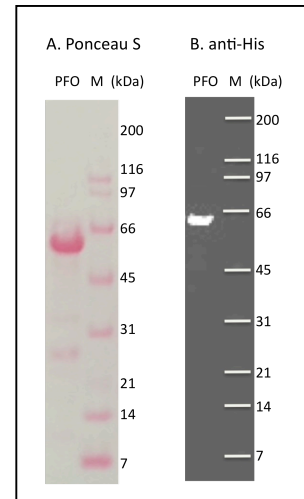
Endotoxin Untested

Formulation 20mM HEPES-Na (pH7.9), 100mM NaCl, and 1mM EDTA.

Storage The protein sample can be stored under sterile conditions at 2- 8°C for one month or at -70°C for three months without significant degradation.

Avoid repeated freeze-thaw cycles.

FOR RESEARCH USE ONLY!!!



Recombinant 6H-PFO* (Y181A,C459A)

Appendix I: Validated seq of 6H-PFO in pET28a vector

```
tttgtttaactttaagaaggagatatacatatgCGGGGTTTctcatcatcatcatcat
      M R G S H H H H H H
ggtatggctagcatgactggtggacagcaaatgggtcgggatctgtacgacgatgacgat
G M A S M T G G Q Q M G R D L Y D D D D
aaggatccgagctcgagatccaaggatataacagataaaaatcaaagtattgattctgga
K D P S S R S K D I T D K N Q S I D S G
atatcaagtttaagttacaatagaaatgaagtttttagctagtaatggagataaaattgaa
I S S L S Y N R N E V L A S N G D K I E
agctttgttccaaaggaaggtaaaaaaactggtaataaatttatagttgtagaacgtcaa
S F V P K E G K K T G N K F I V V E R Q
aaaagatcccttacaacatcaccagtagatatatcaataattgattctgtaaattgaccgt
K R S L T T S P V D I S I I D S V N D R
acatatccaggagcattacaacttgcagataaagcatttgggaaaatagacctacaatc
T Y P G A L Q L A D K A F V E N R P T I
ttaatggtaaaaagaaagcctattaacattaatatagatttaccaggattaaagggcgaa
L M V K R K P I N I N I D L P G L K G E
aatagtataaaggttgatgatccaacctatggaaaagtttctggagcaattgatgaatta
N S I K V D D P T Y G K V S G A I D E L
gtatctaagtggaatgaaaagtattcatctacacatactttaccagcaagaactcaagct
V S K W N E K Y S S T H T L P A R T Q A
tcagaatctatggtttatagtaaatacacaatatcaagtgcccttaattgtgaatgctaaa
S E S M V Y S K S Q I S S A L N V N A K
gtccttgaaaactcacttggagtagactttaatgcagtagcaaacaatgagaaaaaagtt
V L E N S L G V D F N A V A N N E K K V
atgatttttagcatataaacaatatctatacagtaagtgcagatttacctaagaatcca
M I L A Y K Q I F Y T V S A D L P K N P
tcagatctttttgatgacagtgttacatttaattgatttaaaacaaaaggaggtaagtaat
S D L F D D S V T F N D L K Q K G V S N
gaagcacctccacttatggtttcaaagttagcttatggaagaacaatatatgttaagtta
E A P P L M V S N V A Y G R T I Y V K L
gaaactacttctagtagtaaaagatgtacaagctgctttcaaagctcttataaagaact
E T T S S S K D V Q A A F K A L I K N T
gatataaaaaatagtaacaatataaagatatttatgaaaatagttccttcacagcagta
D I K N S Q Q Y K D I Y E N S S F T A V
gttttagaggagatgcacaagaacataacaaagttgtaactaaagacttttgatgaata
V L G G D A Q E H N K V V T K D F D E I
agaaaagtaattaaagacaatgcaacttttagtacaaaaaacccagcatatccaatatct
R K V I K D N A T F S T K N P A Y P I S
tatactagtgttttcttaaaagataactcagttgctgctgttcacaataaaacagattat
Y T S V F L K D N S V A A V H N K T D Y
atagaacaacttctacagagtattctaagggaaaaataaacttagatcatagtggagcc
I E T T S T E Y S K G K I N L D H S G A
tatgttgcacagtttgaagtagcctgggatgaagtttcatatgacaaagaaggaaatgaa
Y V A Q F E V A W D E V S Y D K E G N E
gttttaactcataaaacatgggatggaattatcaagataaaacagctcactattcaaca
V L T H K T W D G N Y Q D K T A H Y S T
```

Recombinant 6H-PFO* (Y181A, C459A)

gtaatacctcttgaagctaatgcaagaaatataagaataaaaagcaagagaggctacaggc
V I P L E A N A R N I R I K A R E A T G
cttgcttgggaatggtggagagatggtataagtgaatatgatggtccattaacaaataat
L A W E W W R D V I S E Y D V P L T N N
ataaatgtttcaatatggggaacaactttataccctggatctagtattacttacaattaa
I N V S I W G T T L Y P G S S I T Y N *
gaattc