

PRMT8 T (Protein arginine N-methyltransferase 8; HRMT1L3, HRMT1L4)

R
R
C
T
1

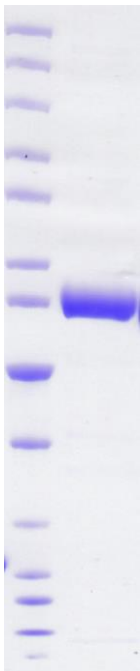
S R P T N Human recombinant PRMT8 (residues 1-344 (N-terminus); C-terminal A fusion protein with N-terminal ST fusion protein and C-terminal Strep-tag, in *Sf9 insect cells*. *9* *a. P* *a t e* *ar inine*
meth ltransferase catal es the transfer of a meth l rou from S a enos I L methionine S to an ω

residues 1 ^ω dis i l t i e i t t e s u s t r t e T R. . RBC's PRMT8 is deleted for

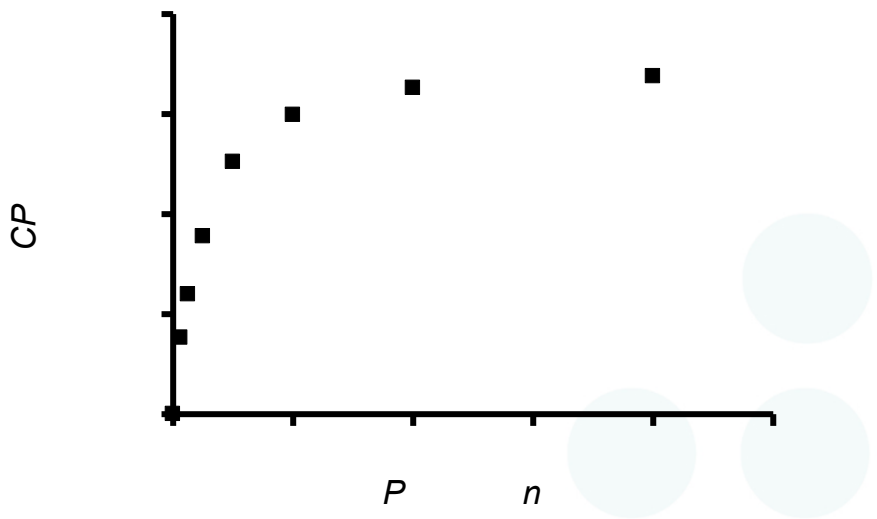
RBC's PRMT8 dis l s et l t r s f e r s e t i t t e e o e t r t i o s o f 1 . M d o e i t T
 R T fused to fill residues 1-8 d e o s l e t i o i e s s u s t r t e s . t i t s d e t e r i e d s T C
 re i t t e d o u t s i s i t t i o f i l t e r l t e s s M u l t i s r e e B T o u t . R e t i o o d i t i o s M T r i s C l 8 . M
 C l M M C l 1 M T T C i . i t s u s t r t e s s i d i t e d o e .

P P .1 l i M T r i s C l 8 M C l 1 l e r o l . M T C P M T T s d e t e r i e d 8
 T R C . T u i l d s t o r e o i e e f o r e u s e . T e r e i i u s e d u d i l u t e e s o u l d e r e f r o e u i l
 f o r e l e s f r e e i i d r i e t o l t o r l i u i d i t r o e . r e e i d s t o r e o f d i l u t e e e i s o t
 r e o e d e d .

R R C 1 . e e e t a l . J . B i o l . C h e m . 9 . . e n e t a l . P L O S O n e e . o u s a a e t a l . N e u r o s c i e n c e 9
 J . S a e h e t a l . J . B i o l . C h e m .



Coomassie blue stained SDS PAGE of PRMT8 T. Markers left are from...



Methylation of P... eth lation determine as C...
 r e c i t a l e c o u n t s i n a s c i n t i l l a t i o n f i l t e r l a t e a s s a . e a c t i o n s e r e
 L m i n . C i t h S a n S a s
 s u s t r a t e s .

This product is not intended for therapeutic or diagnostic use in animals or in humans.