

E. coli Protein Purification

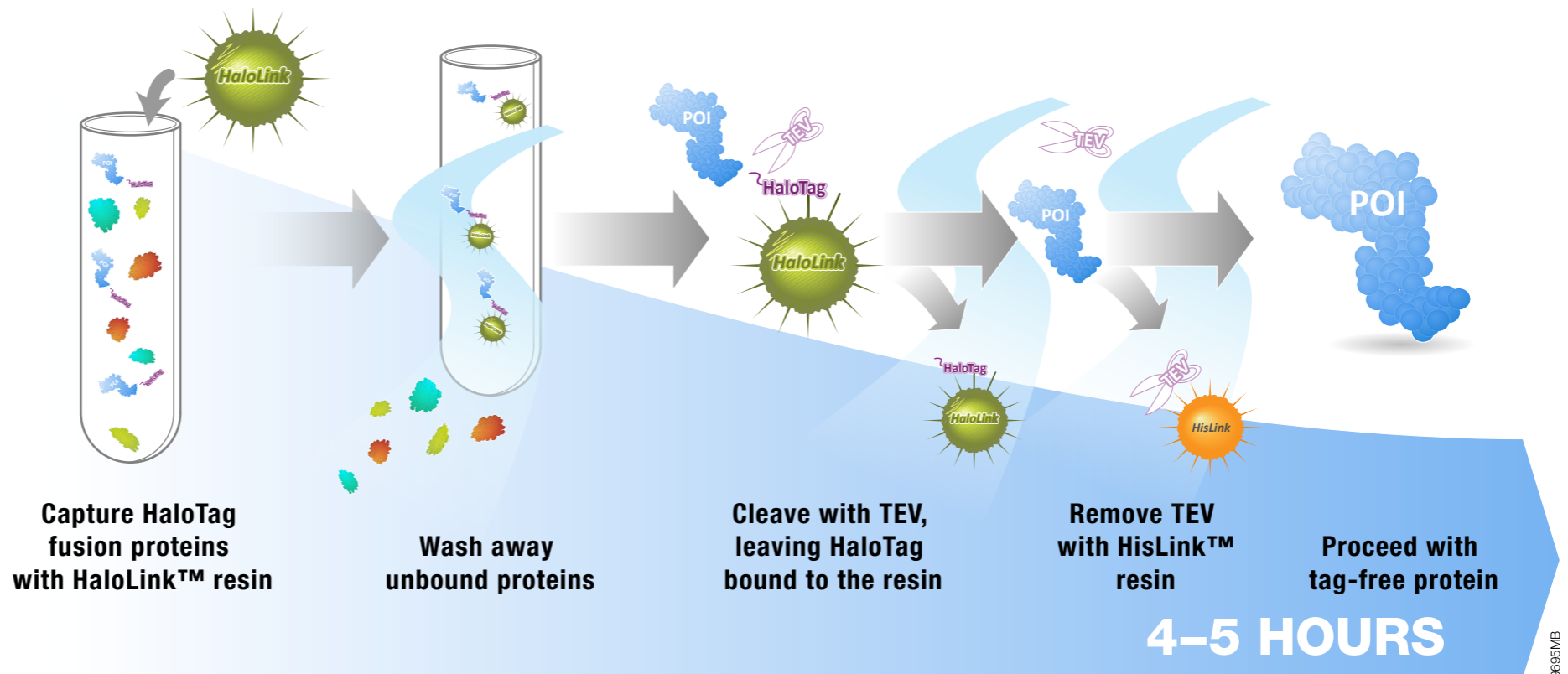
with the HaloTag[®] platform



- ✓ *HaloTag is a highly-effective solubility tag*
- ✓ *Yields more active protein*
- ✓ *Improved protein recovery due to highly specific, covalent binding*

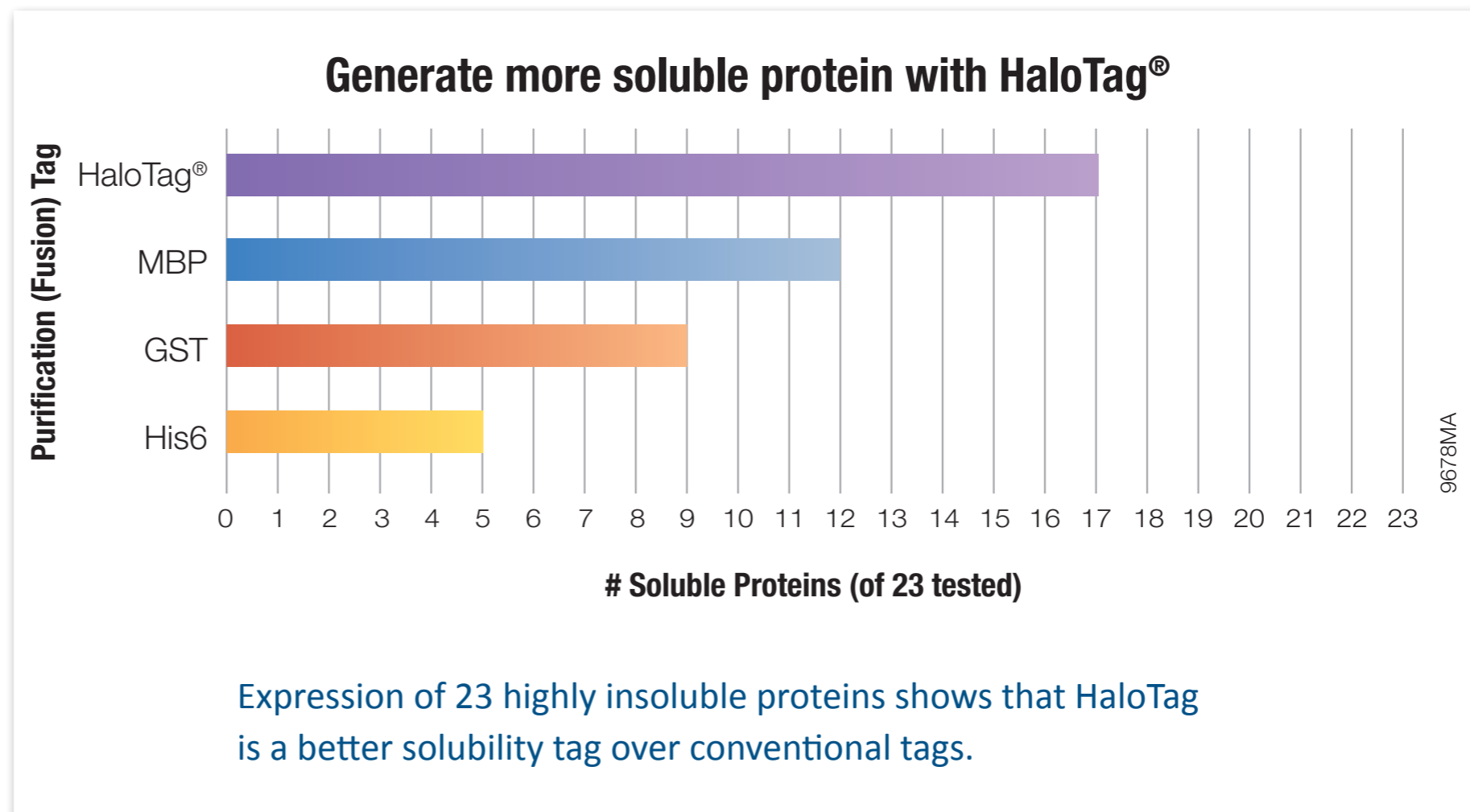


Rapid HaloTag® protocol



This protocol can be further simplified by using an optimized on-column TEV, called HaloTEV protease.

Enhance protein solubility



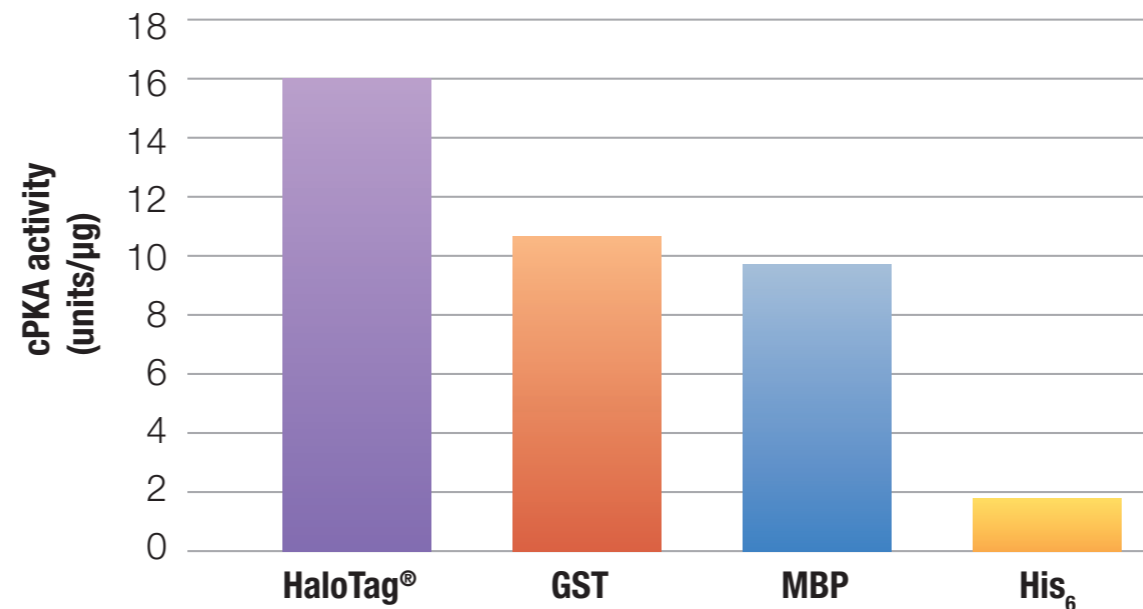
Ohana, R.F. *et al.* (2009) HaloTag7: a genetically engineered tag that enhances bacterial expression of soluble proteins and improves protein purification. *Protein Expr. Purif.* **68** (1): 110-20.

Recover more active protein

HaloTag® purified proteins are highly functional

Gentle purification conditions using a single native buffer and fewer steps in the protocol results in highly active protein

No dialysis necessary!

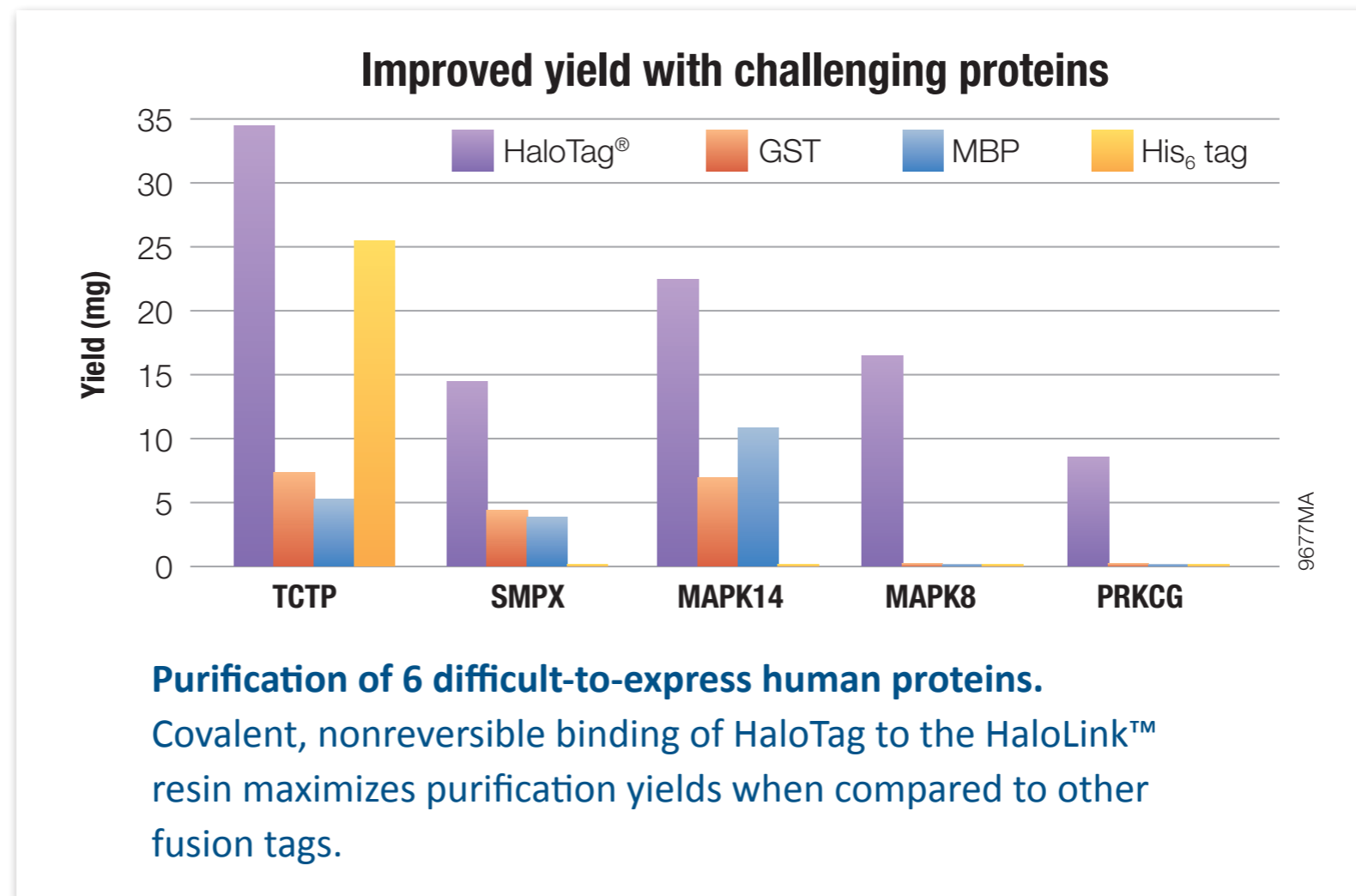


cPKA activity was determined using Kinase-Glo® assay after purification with multiple fusion tags, including HaloTag.

9717MA

Recovery where other tags fail

Higher expression and greater recovery



PROMEGA iSLIDES

FEATURES & BENEFITS

PROTOCOL

EXPRESS & PURIFY
DIFFICULT PROTEINS

PROTEIN ACTIVITY

BETTER RECOVERY

MORE INFORMATION

E. coli PROTEIN PURIFICATION WITH THE HALOTAG® PLATFORM

References

Ohana, R.F. *et al.* (2009) HaloTag7: a genetically engineered tag that enhances bacterial expression of soluble proteins and improves protein purification. *Protein Expr. Purif.*; **68**(1); 110-20.

Hata, T. and Nakayama, M. (2007) Rapid single-tube method for small-scale affinity purification of polyclonal antibodies using HaloTag Technology. *J Biochem. Biophys. Methods*; **70**(4); 679-82.

Motejadded, H. *et al.* (2010) Expression, one-step purification, and immobilization of HaloTag™ fusion proteins on chloroalkane-functionalized magnetic beads. *Appl. Biochem. Biotechnol.*; **162**(7); 2098-110.



Ordering Information

PRODUCT	SIZE	CATALOG #
HaloTag® Protein Purification System	1 each	G6280
HaloTag® Protein Purification System Sample Pack	1 each	G6270

HaloTag and Kinase-Glo are registered trademarks of Promega Corporation. HaloLink is a trademark of Promega Corporation.

