



# Important Notice

for the use of ProteasMAX™ Surfactant, Trypsin Enhancer  
V2071, V2072

Dear customer,

thank you for choosing ProteasMAX™ Surfactant, Trypsin Enhancer to support your trypsin digestion. To facilitate your daily work with ProteasMAX™ please consider the notes we have put together below.

1. ProteasMAX™ Surfactant, Trypsin Enhancer is designed to degrade over the course of a trypsin digestion reaction. We recommend 0.025% ProteasMAX™ concentration for protein digestion. A maximum concentration of 0.05% can be used for in-solution digestion of membrane proteins. Higher concentrations may cause peptide precipitation.
2. 0.2% or even higher ProteasMAX™ concentration can be required to solubilize difficult proteins such as membrane proteins. Please dilute the solubilized proteins to acceptable ProteasMAX™ concentration (see 1.) before digestion.
3. ProteasMAX™ is currently recommended for in-gel digestion of proteins present at amounts detectable with Coomassie Blue R250 (50 ng and up). Application of ProteasMAX™ for in-gel digestion of lower protein amounts is under development.
4. We recommend 0.6 µl C18 ZipTip® or comparable C18 tips (i.e. 10 µl OMIX® Tips) to clean up and concentrate peptides after in-gel digestion. The use of 0.2 µl C18 tips can lead to a significant peptide loss due to competition of the degraded surfactant and peptides for binding sites.
5. At sufficiently high protein amounts the digested sample can be directly analyzed with MALDI without SPE. ProteasMAX™ degradation products do not interfere with MALDI.
6. Direct analysis of the digested sample by ESI is not recommended. Analysis by ESI should be preceded by LC or SPE.

For further details please carefully read the Technical Bulletin 373 which accompanies this product shipment. Should you wish personal assistance please contact Dr. Jan Adam from Technical Service at 0621-8501-278.

We wish you a successful MS sample preparation!  
Sincerely,  
Promega GmbH