

MY FAVORITE PROMEGA PRODUCT TESTIMONIAL I Like, I Share!



Promega remercie tous les participants de leurs témoignages lors du dernier congrès EuPa 2013 !

Que 2014 soit l'occasion de toujours plus d'échanges et de rencontres riches en expériences !

Trypsin Gold *"Very reproductible digestion!"*

Magalie, Institut Pasteur

Trypsin/lys C mix *"Very satisfied"*

François, SANOFI

Trypsin *"Efficient & reproductible, easy to use, I like!"*

Edith, DCMR Ecole Polytechnique*

Trypsin Gold *"The best enzyme I know for a short time of digestion (03:00)!"*

Géraldine, CLIPP Proteomic Platform

Trypsin Gold *"Never had any bad surprise with this high quality trypsin"*

Virginie, LEBS

Trypsin *"Great analysis for imagerie in clinical proteomic field"*

Claude, PITZ CRUZ Faculté de Pharmacie

Trypsin, AspN, ProteaseMax *"We use to process our protein samples for protein identification via trypsin, together sometimes with Proteasemax if membranes proteins. We also use AspN for completing peptide mass fingerprint"*

Régine, Proteomic Platform

Trypsin Gold *"As an engineer on a proteomic platform, I am very satisfied with the quality of this product routinely used for our protein digestion (in gel or in solution)"*

Martial, IGF

Trypsin sequencing grade *"For many years now, the proteomic core facility biogenouest purchases several promega products because of the high quality of their products fit perfectly with our process and tehcnology platform to provide superior results"*

Régis, BIOGENOUEST

Western blot ECL *"Western blot in the west of France with Promega!!!"*

Charlotte, IFREMER

ProteaseMax *"It's a good product for in gel and in solution trypsin digestion"*

Valérie, INRA

ProteaseMax *"Improving sequence coverage in proteomic is a must"*

Julien, INRA

Trypsin *"The best enzyme for protein digestion until the beginning of my proteomics experiments 12 years ago"* Emmanuelle, BIOGENOUEST

Trypsin *"Good product"* Samuel, IBPC

Trypsin *"Always satisfied about quality of Promega trypsin. Efficiency"*

Jessica, PISSARO Université Rouen

Trypsin *"Very reproducible"*

Catherine, SANOFI

Trypsin sequencing grade *"Very good reproductibility intra batch"*

Armelle, SANOFI

Trypsin *"Great analysis for imagerie in clinical proteomic field"*

Claude, PIZ CRUZ Faculté de Pharmacie

Trypsin Gold *"Great product reproductibility"*

Sylvie, IBP Plateform PROMETHEE

Trypsin sequencing grade *"Your MS specific modified trypsin is much better than other trypsin fo in solution digest : less autolysis"*

Philippa, INRA

Trypsin *"Good results reproductibility efficiency"*

Mathilde, IGBMC

Trypsin Gold *"Great easy to use and effective, thanks!"*

Marven, CLIPP Proteomic Platform

Trypsin Gold *"Good reproductibility and good conservation"*

Aurélie, INRA

Trypsin *"Excellent product reproductible"*

Stéphane, CRCM UMR 1068

Trypsin *"Easy to use, reproductible results"*

Sophie, DCMR Ecole Polytechnique

Trypsin *"Efficient"*

Marie, MNHN

Trypsin *"I always use trypsin for my protein digestion"*

Mathilde, INRA

Trypsin Gold *"Great enzyme for proteomic analysis"*

Emmanuelle, IPBS

Trypsin *"Works very well for digestion of protein before identification or MALDI mapping of peptides"*

Mélanie, BIOGENOUEST

Gotaq qPCR master mix *"Good product, good price"*

Laetitia, INRA

Gotaq qPCR master mix *"Good product"*

Aurélie, INRA