

Offered in Europe

Pre-paid

1 Sample

5 Samples

91-10x

Two specific part-codes are available  
depending on the quantity of  
samples to be tested.

95-10x

91-101

Typically water  
sample

Tested 1/1  
by default

91-102

Typically dialysis  
sample

Tested 1/5 and  
1/10 by default

91-103

Typically research  
sample

Tested 1/10,  
1/100 and 1/1000  
by default

Dilution scheme

Three options are  
available depending  
on the quantity of  
dilutions to be tested

95-101

Typically water  
sample

Tested 1/1  
by default

95-102

Typically dialysis  
sample

Tested 1/5 and  
1/10 by default

95-103

Typically research  
sample

Tested 1/10,  
1/100 and 1/1000  
by default

80-508

Special treatment

The Pre-paid Services do not include special handling on your product. On request, a specific treatment is applied or special reagents are used (pH adjustment, heat inactivation, use of Tris-Buffer,  $MgCl_2$ , Pyrosperser™ or glucan blocker.)