

## HistaStrip<sup>™</sup> Test Kits

CATALOG #FOOD-1100-01 (24 Tests)

The HistaStrip™ Test Kit is a convenient dipstick technology used for the rapid visual detection of histamine in seafood, fish meal, milk, and fish sauce. The test strips can detect unsafe levels of histamine in only 4 minutes without the use of organic solvents or heating steps.

The HistaStrip™ test strips contain a pad embedded with enzyme that reacts specifically with histamine to chemically reduce a dye indicator, producing a color change for easy visual interpretation of results. When the strips are dipped into samples containing histamine, the pad rapidly changes from white to a color designating the histamine concentration present in the food or drink. Detection capabilities using the strips are well below global Maximum Residue Limits (MRL). The self-contained, visual test enables easy testing in and out of the laboratory.

The HistaStrip<sup>™</sup> can also give full quantitative results using the QuickSTAR<sup>™</sup> Strip Reader, enabling the customer to determine exact levels of histamine in the sample and easily save data to a computer and quality assurance system.

## The features of the kit are:

- 4 Minute results
- Visual semi-quantitative test
- Full quantitative results with reader

Sample:					
Seafood	0	25	50	75	100
Milk	0	12	24	36	48
Fish Sauce	0	125	250	375	500
Fishmeal	0	500	1000	1500	2000

PerkinElmer, Inc. 7050 Burleson Road Austin, TX 78744 USA P: (888) 208-2246 or (+1) 512-707-8993 www.biooscientific.com



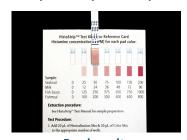
Homogenize sample



Add enrichment solution & mix



Dip test strip into sample



**Read results** 

