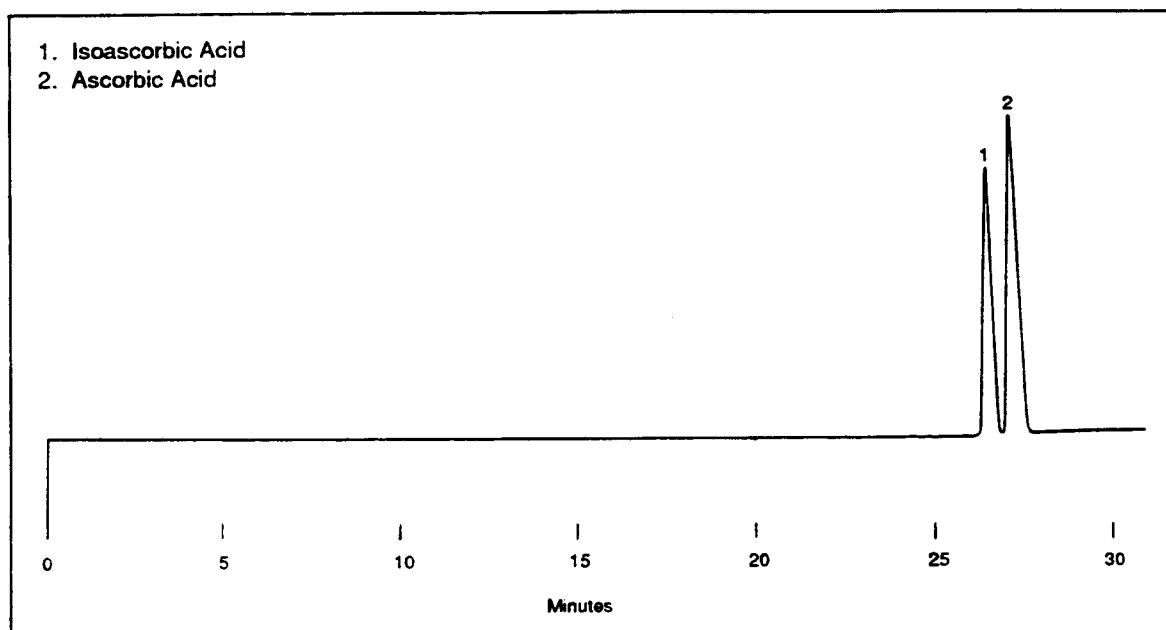


Ascorbic Acid (Vitamin C) and Isoascorbic Acid (Erythorbic Acid)Application
Note: 15

Ascorbic acid is a water-soluble vitamin with a wide natural distribution. Isoascorbic acid, its stereoisomer, has only 5% of the biological activity of ascorbic acid. Because of its similar redox potential, it is commonly used in preference to ascorbic acid as an antioxidant in foods. Capillary electrophoresis can be used to separate these two compounds.



Capillary:	50 cm x 50 μ m, coated
Buffer:	0.1 M sodium phosphate, pH 5.7
Load conditions:	8 kV, 5 seconds
Run conditions:	8 kV, constant voltage, $\ominus \rightarrow \oplus$ polarity
Detection:	UV, 265 nm, 0.02 AUFS