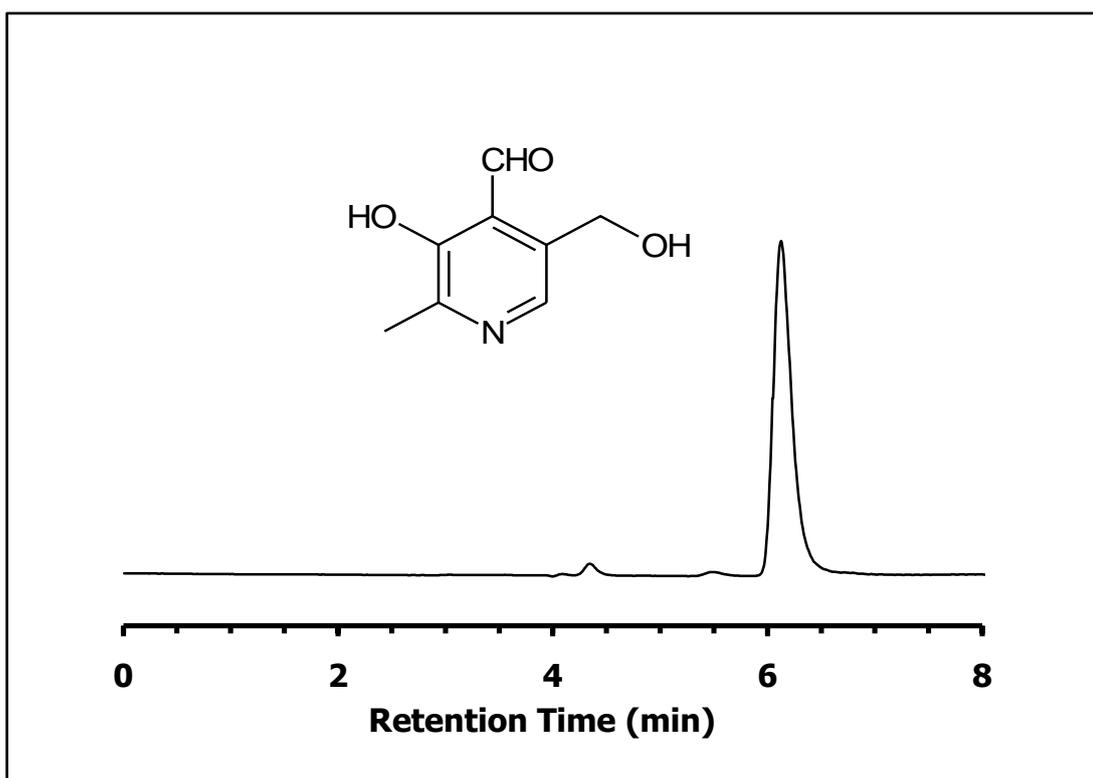


ZIC®-HILIC Separation of Pyridoxal

Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 150 x 4.6 mm, 5 µm, 200 Å (cat. no. 150455)
Injection: 10 µL in mobile phase
Detection: UV @ 254 nm (UFS 1.0 V)
Pressure Drop: 1.6 MPa (230 psi)
Flow Rate: 0.5 mL/min
Mobile Phase (v/v): 90%, Acetonitrile
10%, Ammonium Formate 10 mM
(total ionic strength = 1 mM)



Chromatographic Data

| No. | Compound | | Time (min) | Retention factor |
|-----|------------|-------------|------------|------------------|
| | t_0 | void volume | 3,0 | - |
| 1 | Vitamin B6 | Pyridoxal | 6,1 | 1,0 |