



- | Analytes Included | Quantify Up to 11 Routinely Tested Cannabinoids | | | | | | | Differentiate Acids & Neutrals | | | |
|---------------------|---|---|--|---|---|---|---|---|---|---|---|
| | 11-Mix | 10-Mix | 9-Mix (CA) | 6-Mix | 5-Mix | 4-Mix | 3-Mix | Neutrals 8-Mix | Neutrals 9-Mix | Acids 6-Mix | Acids 7-Mix |
| CBD | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | | |
| Δ ⁹ -THC | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | | |
| CBN | ✓ | ✓ | ✓ | ✓ | ✓ | | ✓ | ✓ | ✓ | | |
| THCA-A | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ | | | | ✓ | ✓ |
| CBDA | ✓ | ✓ | ✓ | ✓ | ✓ | | | | | ✓ | ✓ |
| CBG | ✓ | ✓ | ✓ | ✓ | | | | ✓ | ✓ | | |
| CBC | ✓ | ✓ | ✓ | | | | | ✓ | ✓ | | |
| Δ ⁸ -THC | ✓ | ✓ | ✓ | | | ✓ | | ✓ | ✓ | | |
| CBDV | ✓ | ✓ | | | | | | ✓ | ✓ | | |
| THCV | ✓ | | ✓ | | | | | ✓ | ✓ | | |
| CBGA | ✓ | ✓ | | | | | | | | ✓ | ✓ |
| CBDVA | | | | | | | | | | ✓ | ✓ |
| THCVA | | | | | | | | | | ✓ | ✓ |
| CBCA | | | | | | | | | | ✓ | ✓ |
| CBL | | | | | | | | | ✓ | | |
| CBNA | | | | | | | | | | | ✓ |
| Mixture Diluent | acetonitrile | | | | | | methanol | acetonitrile | methanol | 99:1 acetonitrile:DIPEA | |
| Fill Volume | 0.5 ml or 1 ml | 1 ml | | | | | | 0.5 ml | 1 ml | 0.5 ml | 1 ml |
| Available Formats | View Online  | View Online  | View Online  | View Online  | View Online  | View Online  | View Online  | View Online  | View Online  | View Online  | View Online  |