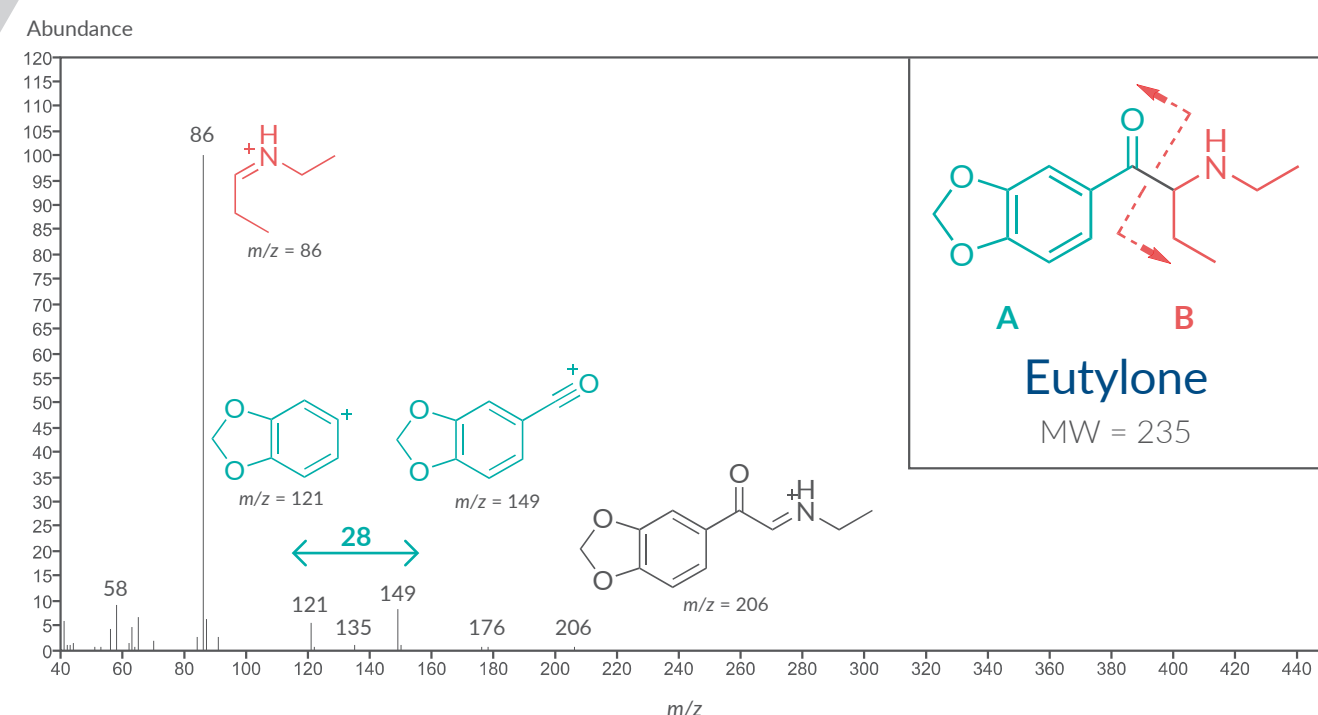
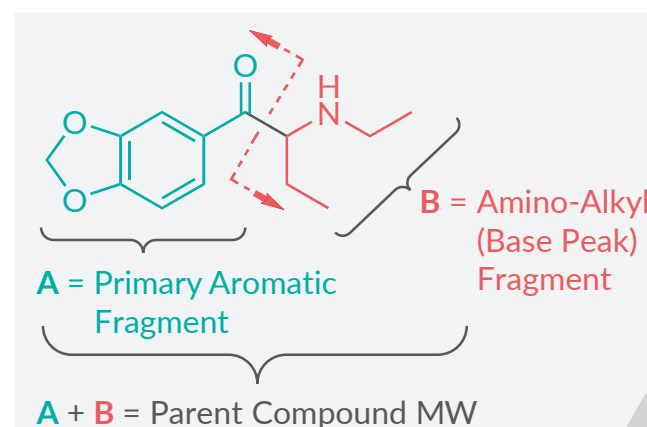


Mass Spectrum of Eutylone and Tips for Interpretation



Tips for EI-MS Interpretation:

- Fragmentation often results in no molecular weight ion.
- Fragmentation occurs via α -cleavage to yield the **primary aromatic (A)** and **amino-alkyl (B)** fragments.
- The sum of the **A** and **B** fragments results in the parent compound molecular weight (MW).
- Base peak is always the **amino-alkyl fragment (B)**, which may also be observed in amphetamines.
- Fragment A**, the acylium ion, further cleaves to yield a subsequent fragment minus 28 (**A-28**), also common to amphetamines.
- IR spectroscopy and/or NMR can be used in addition to GC-MS to differentiate amphetamines from cathinones.
- Multiple positional isomers are possible for cathinones.



Laboratory Guide for CATHINONE Identification and Naming

Common Substitutions

PHENYL RING

R_1 denotes hydrogen or one or more substitutions on the phenyl ring (alkoxy groups, halides, alkyl/aryl groups, etc.)

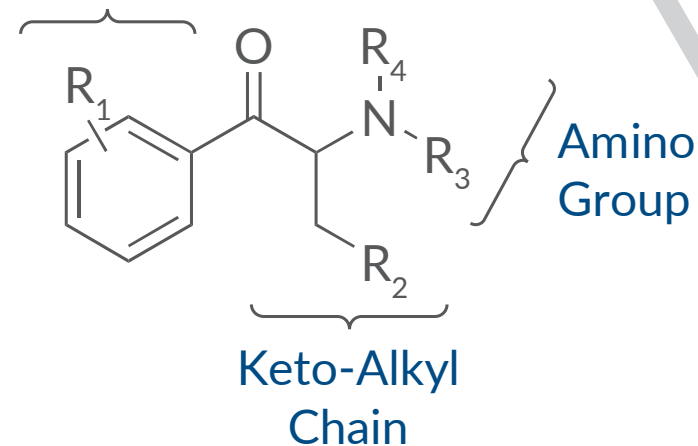
KETO-ALKYL CHAIN

R_2 denotes hydrogen or increased carbon chain length, which can also be branched

AMINO GROUP

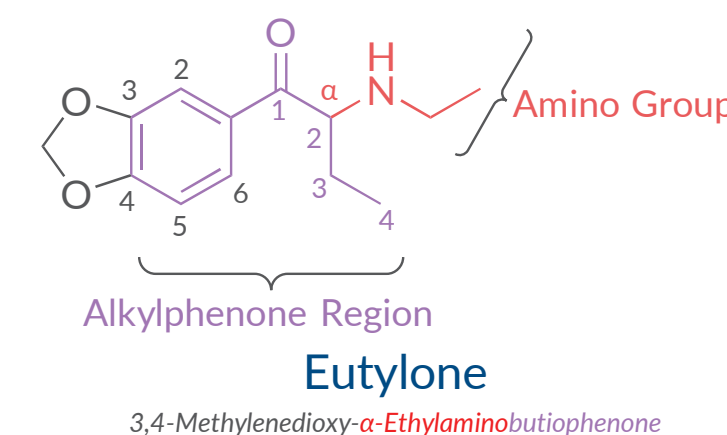
R_3/R_4 denotes hydrogen, alkyl groups, or a ring system

Phenyl Ring



Standardized Alkylphenone Naming System

- The first step in using this naming system is recognition of the **alkylphenone region**. Alkylphenones are named "propiofenone," "butiofenone," "pentio/valerophenone," etc. based on the keto-alkyl chain length (see [Table](#)).
- The **α -amino group** is then identified, such as " α -methylamino," " α -ethylamino," etc. The α -amino group name is positioned before the alkylphenone name in the naming order sequence. If the amine group is a ring system such as a pyrrolidine or piperidine, then " α -pyrrolidino" and " α -piperidino" are used.
- Substituents on the phenyl ring and their location designation ("4-fluoro," "3,4-methylenedioxy," etc.) are positioned at the beginning of the name.



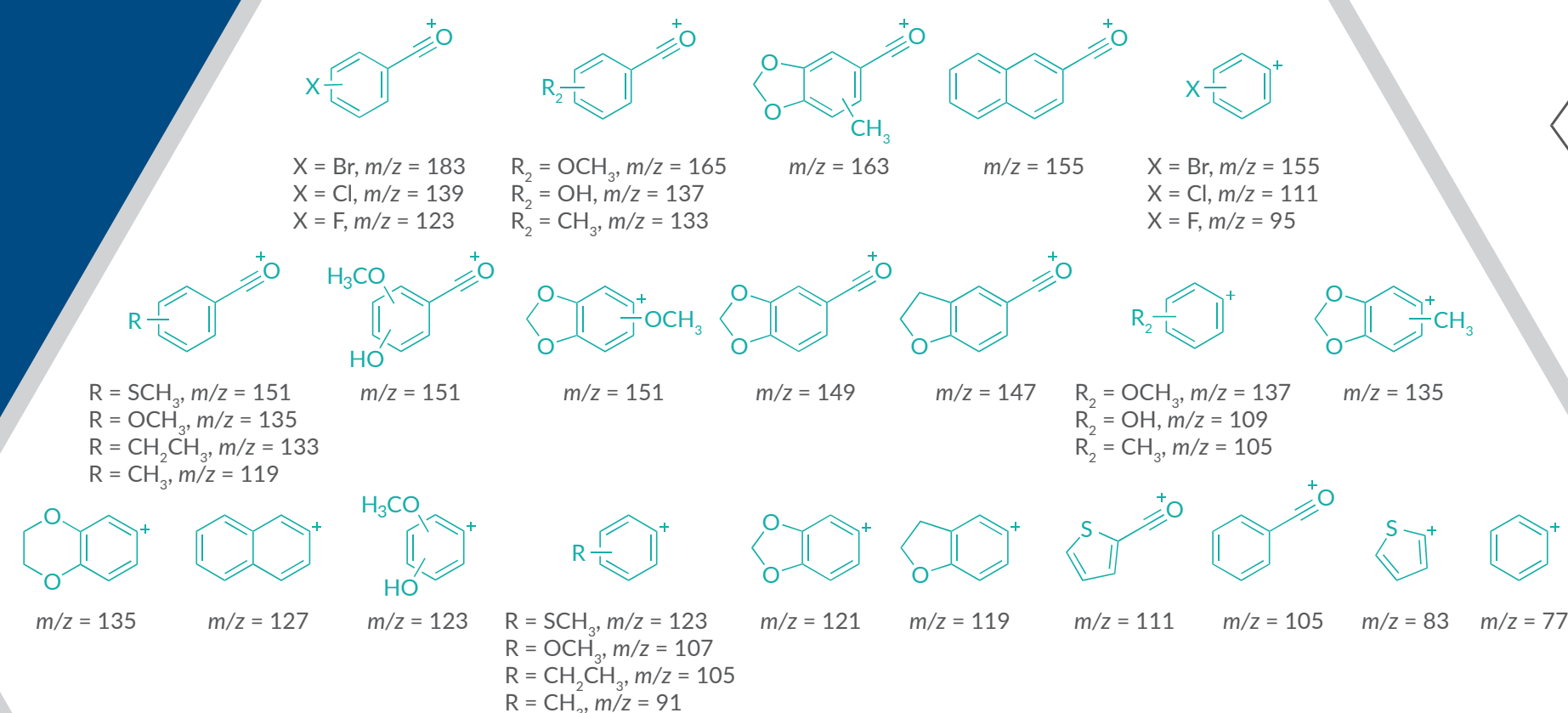
Alkylphenone Region Naming	Keto-Alkyl Chain (# of Carbons)
Propiofenone	3
Butyro/Butano/Butiofenone	4
Pentio/Valerophenone	5
Hexio/Hexanophenone	6
Heptanophenone	7
Octanophenone	8
Nonanophenone	9
Decanophenone	10

Final Naming Order Sequence: Phenyl ring substituents then **α -amino group** then **alkylphenone region**

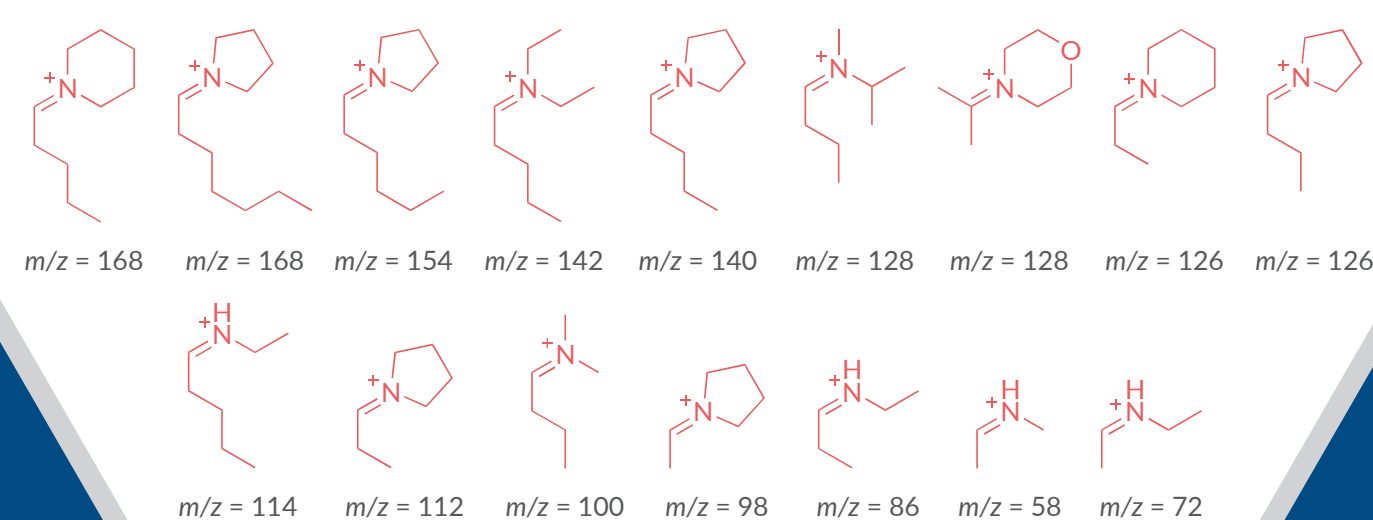
Examples: 3-fluoro- α -Ethylaminobutiofenone or 4-methyl- α -Isopropylaminovalerophenone

Common Mass Spec Fragments

Aromatic Fragments (in descending m/z order)

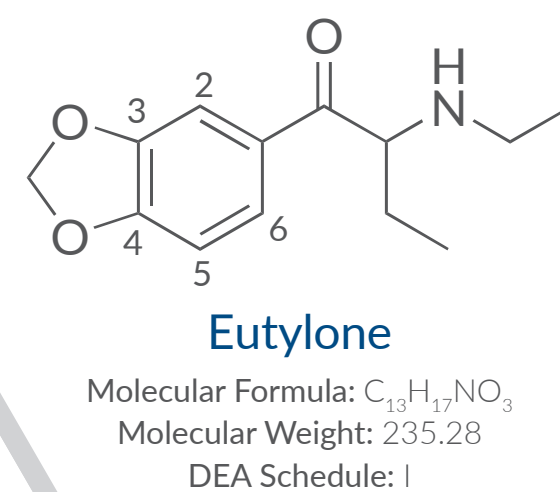


Amino-Alkyl (Base Peak) Fragments (in descending m/z order)



Please note that multiple isomers can exist for all mass spec fragments.

Substituted Cathinone Example: Eutylone



SYNONYMS

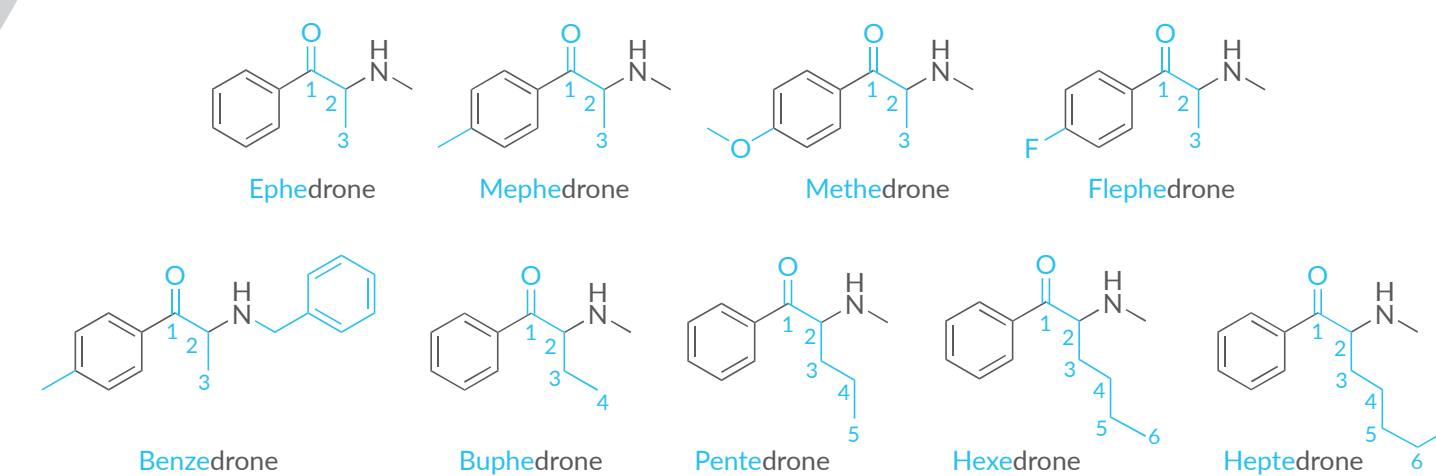
β -keto-Ethylbenzodioxolylbutanamine,
3,4-Methylenedioxy- α -Ethylaminobutiofenone,
bk-EBDB, "Ethyl Butylone"

SUBSTITUTIONS

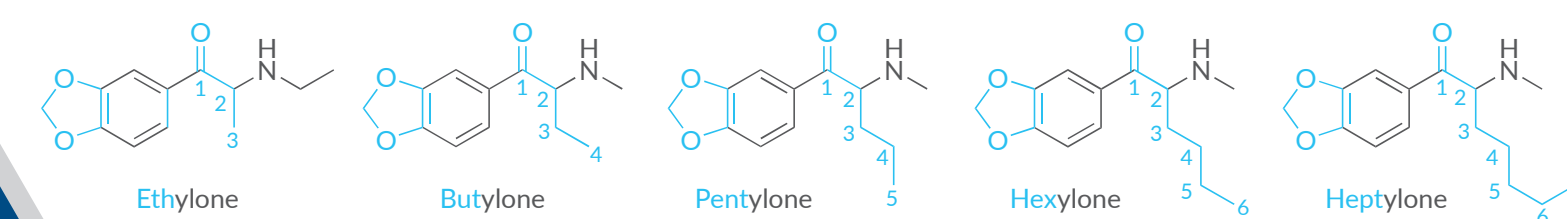
Contains a methylenedioxy on the phenyl ring; R_2 is a methyl group, R_3 is an ethyl group, and R_4 is a hydrogen

Other Common Naming Conventions

"drone" Naming: The suffix "drone" commonly represents a cathinone without a methylenedioxy group, and the prefix typically represents the length of the alkyl chain. Please note that there are exceptions to this trend such as in the case of the first five examples below.



"ylone" Naming: The suffix "ylone" commonly represents a cathinone with a methylenedioxy group fused to the aryl ring, and the prefix typically corresponds to the length of the alkyl chain.



β -keto Naming: Some cathinones are named after their corresponding amphetamine analogs as " β -keto" (bk) versions.

