

# Benzo[a]pyrene

Polycyclic aromatic hydrocarbon (PAH)  
found in coal tar, cigarette smoke, and wood  
smoke; carcinogen.

Environmental carcinogen found in automobile exhaust, jet engine exhaust, coal tar, tobacco smoke, etc. It is carcinogenic to laboratory animals.

## Ordering Information

[Order Online »](#)

LKT-B1652-M500	500mg
LKT-B1652-G001	1g

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Alternative Name**      1,2-Benzpyrene, 1,2-Benzopyrene, 6,7-Benzopyrene, 3,4-Benzylpyrene, 3,4-Benz[a]pyrene

**CAS**      50-32-8

**Formula**       $C_{20}H_{12}$

**MI**      **13:** 1104

**MW**      252.32

**Purity**       $\geq 98\%$

**Solubility**      Soluble in chloroform, benzene, toluene, xylene, ether. Insoluble in water. Sparingly soluble in alcohol, methanol.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)