

Bourgeonal

hOR17-4 ligand

A previously undescribed human testicular olfactory receptor known as hOR17-4 was cloned and functionally expressed. Bourgeonal is a potent agonist at hOR17-4 and acts as a strong chemoattractant in sperm behavioral assays. The bourgeonal, hOR17-4 signaling pathway potentially governs chemical communication between sperm and egg and may be used to manipulate or block fertilization.

Citations: 3

[View Online »](#)

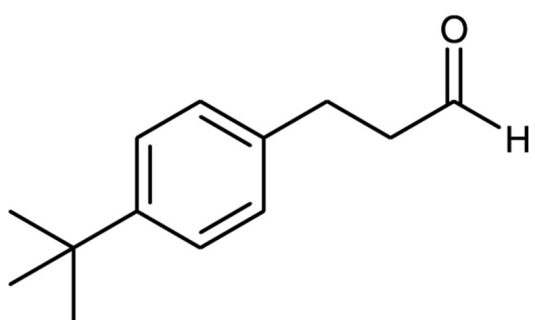
Ordering Information

[Order Online »](#)

BML-N156-0025	25mg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-t-Butylbenzenepropionaldehyde
Appearance	Colorless oil.
CAS	18127-01-0
Couple Target	hOR17-4
Couple Type	Ligand
Formula	$C_{13}H_{18}O$
MW	190.3
Purity	≥90% (TLC)
Solubility	Soluble in DMSO (25 mg/ml) or ethanol (25 mg/ml).



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)