

6-CP-FAD

Functionalized analogue of FAD

6-CP-FAD is a functionalized analogue of the adenine-containing enzymatic redox cofactor flavin adenine dinucleotide (FAD). The terminal carboxy group of 6-CP-FAD can be used for coupling of various labels, dyes or supports with free amino groups. Via its corresponding succinimidyl ester, 6-CP-FAD can also be coupled to FAD-binding apo-proteins.

BLG-F017-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

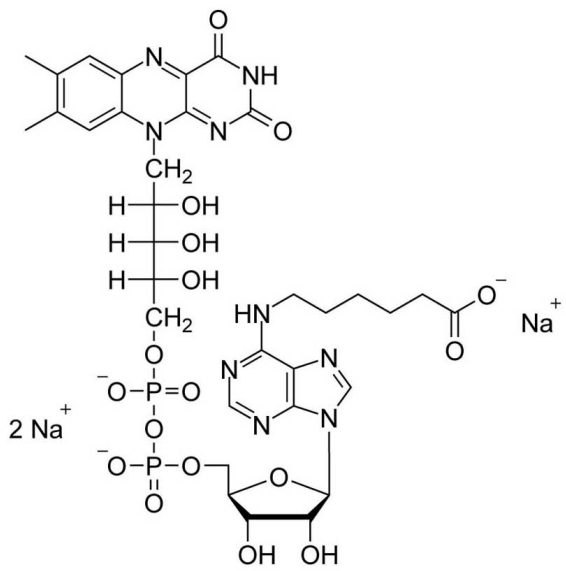
Ordering Information

[Order Online »](#)

BLG-F017-01	1 μ mol
BLG-F017-05	5x1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling Protect from light.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Flavin-N6-(5-carboxypentyl)adenine dinucleotide . sodium salt

CAS 177718-13-7

Couple Type Modified nucleotides

Formula $C_{33}H_{43}N_9O_{17}P_2$ (free acid)

MW 899.7 (free acid)

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.9 mg



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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